

Testing can also be ordered via online portal – please scan or click on QR code.  
Please complete every field and tick box clearly.



STEP 1: PATIENT INFORMATION

Patient's First Name Middle Initial Patient's Last Name  
     
 Patient's Date of Birth Patient ID/MR Number/External Sample Number  
 Biological Sex:  Male  Female  Unknown  
 Gender Identity (if different from above):   
   
 Patient's Street Address City / Town  
      
 State Zip Code Country Patient's Preferred Phone Patient's Email  
 Ethnicity (check all that apply):  African-American  Asian (China, Japan, Korea)  Caucasian/N. European/S. European  Finnish  French Canadian  Hispanic  
 Jewish - Ashkenazi  Jewish - Sephardic  Mediterranean  Middle Eastern (Saudi Arabia, Qatar, Iraq, Turkey)  Native American  E. Indian  
 Southeast Asian (Vietnam, Cambodia, Thailand)  South Asian (India, Pakistan)  Other (specify)

PATIENT SAMPLE INFORMATION

**SAMPLE TYPE:**  Whole Blood  Saliva Swab  Urine  Dried Blood Spots  C-Dried Blood Spots  DNA, Source:   DHeparin - Plasma  EDTA - Plasma  Other:   
 Collection Date: MM/DD/YY  
 Was this sample collected in the State of NV, NY or OR?:  Yes  No  
 (If yes, separate consent is required. See forms section of website.)

INDICATION FOR TESTING

Clinical Diagnosis:  (medical records/clinical notes are required.) Age at Initial Presentation:

STEP 2: ORDERING PROVIDER AND REPORTING PREFERENCES

Provider's First and Last Name NPI  
   
 Clinic/Hospital/Institution Name Provider's Email  
      
 Provider's Street Address City / Town State Zip Code Country  
    
 Provider's Phone Provider's Fax How would you like to receive the report?:  
 Fax  Email  Portal

SEND ADDITIONAL COPY OF RESULTS TO (If applicable)

Name Role with patient/Job title Clinic/Hospital/Institution Name  
     
 Phone Number Fax Number Email Address How would you like to receive the report?:  
 Fax  Email  Portal

STEP 3: BILLING INFORMATION

INSTITUTIONAL BILLING

Institution/Organization Name Billing Account ID P.O. Number (if applicable)  
   
 Contact Name Contact Phone

PATIENT (SELF) PAYMENT

By providing payment information, you are authorizing Revvity Omics to process payment at the associated charge for tests ordered. Test cost is available on our website, or may be confirmed by calling 877-475-4436. Payment is required prior to test initiation. The patient's sample will be placed on hold (for up to 30 days) until payment is secured. If the patient does not provide payment to Revvity Omics within 30 days, the test order may be canceled. Please note that failure by the patient to respond in a timely fashion to Revvity Omics attempts to obtain payment may cause a delay in the receipt of the results report.

CREDIT CARD (Please fill out all information below)  CHECK: \$  Amount Enclosed (Please make checks payable to: Revvity Omics, Inc.)  
       
 Credit Card Number Exp. Date CVV Cardholder Printed Name as Appears on Card Amount  
      
 Credit Card Billing Street Address City / Town State Zip Code  
   
 Cardholder Signature Cardholder Phone

CONTACT FOR PAYMENT INFORMATION  
    
 Name Phone Email Address

FOR INTERNAL USE ONLY

Date Rec'd	Rec'd			
TEMP	SPEC	COL	#TUBES	VOL
R/C/F				
R/C/F				
R/C/F				

## General Biochemical and Molecular Requisition Form

### STEP 4: TEST MENU

#### BIOCHEMICAL TESTS

##### SCREENING PANELS

- B0200 StepOne® Comprehensive Biochemical Profile\*
- Birth Time: \_\_\_\_\_ Collection Time: \_\_\_\_\_
- Weeks' Gestation: \_\_\_\_\_ Birth Weight: \_\_\_\_\_
- Transfusion status:  Yes  No
- If yes, transfusion type:  Platelet Date: \_\_\_\_\_ Time: \_\_\_\_\_
- Plasma Date: \_\_\_\_\_ Time: \_\_\_\_\_
- RBC Date: \_\_\_\_\_ Time: \_\_\_\_\_
- B0210 Acylcarnitine Profile
- B2020 Amino Acid Profile
- B2040 Lysosomal Storage Disease Enzyme Panel
- B0024 Post-Mortem Screening Panel
- BG100 Nicotinamide Adenine Dinucleotide (NAD) - C-DBS required\*
- GED1D Mucopolysaccharidosis type I (MPS I) marker - DBS required
- GED2D Mucopolysaccharidosis type II (MPS II) marker - DBS required
- CRIMPW Pompe-CRIM analysis

##### DIAGNOSTIC AND MONITORING PANELS

- B0009 Galactosemia Monitoring
- B0018 PKU Clinical Monitoring
- B0022 Tyrosinemia Monitoring

#### COMPREHENSIVE NEWBORN TESTING

- D3005 NeoSeq Newborn and Pediatric Gene Testing
- D3004 Expanded Newborn Screening (NBS) Gene Sequencing Test

##### CURATED PANELS

- D3200 NeoNGS Panel
- D3200F STAT NeoNGS Panel

##### ADDITIONAL TESTING†

- D3100 AnyGene™ Test: Single Gene Sequencing and Del/Dup Test  
Please submit requested gene for testing at [apps-omics.revvity.com/gene-dashboard](https://apps-omics.revvity.com/gene-dashboard), and include custom gene ID below

*Provide gene or custom panel ID here:* \_\_\_\_\_

Test Code: \_\_\_\_\_

Test Name: \_\_\_\_\_

† Additional testing options including DNA Mutation Screens and Gene Sequencing for individual conditions (or sets of conditions) can be found on pages 4 - 7.

\* DBS Required. Test code BG100 requires collection on C-DBS cards.

### STEP 5: PHYSICIAN CONFIRMATION OF INFORMED CONSENT AND MEDICAL NECESSITY

The undersigned person (or designated representative thereof) certifies that: (a) he/she is a licensed medical professional authorized to order the testing ordered herein; (b) he/she fully complies with all applicable federal, state, and local laws, regulations, and rules, including but not limited to those governing genetic testing, informed consent, and patient consent and authorization requirements for the test(s) ordered; (c) he/she will obtain informed consent of the patient in compliance with all applicable laws and regulations, which shall include, to the extent applicable: (i) a statement of the purpose of the test(s) ordered; (ii) a statement that prior to signing the consent form, the consenting person discussed with the medical practitioner ordering the test the reliability of positive or negative test results and the level of certainty that a positive test result for that disease or condition serves as a predictor of such disease; (iii) a statement that the consenting person was informed about the availability and importance of genetic counseling and provided with written information identifying a genetic counselor or medical geneticist from whom the consenting person might obtain such counseling; (iv) a general description of each disease or condition tested for; and (v) the person or persons to whom the test results may be disclosed; (d) he/she will maintain, as part of the patient's record, documentation of the patient's informed consent and authorization for the test(s) ordered that complies with applicable laws and regulations, and will make such documentation available to Revvity upon request; (e) tests ordered are medically necessary and results may impact medical management for the patient; and (f) the information provided on this Test Requisition Form is complete, true, and accurate to the best of his/her knowledge.

Signature \_\_\_\_\_ Date \_\_\_\_\_

## General Biochemical and Molecular Requisition Form

DETAILED MEDICAL RECORDS, PREVIOUS TEST RESULTS AND FAMILY HISTORY MUST BE ATTACHED FOR ALL CASES.  
CLINICAL INFORMATION IS CRUCIAL FOR ACCURATE INTERPRETATION OF RESULTS.

ADDITIONAL PHENOTYPE / PATIENT HISTORY SECTION (Check all that apply)

Clinical diagnosis: \_\_\_\_\_ Age of manifestation: \_\_\_\_\_

### NEUROLOGY

#### 1. Neurodevelopmental abnormality

- 1.1 Autism
- 1.2 Attention deficit disorder
- 1.3 Global developmental delay
- 1.4 Delayed motor development
- 1.5 Delayed language development
- 1.6 Developmental regression
- 1.7 Intellectual disability

#### 2. Brain imaging

- 2.1 Abnormal myelination
- 2.2 Agenesis of corpus callosum
- 2.3 Brain atrophy
- 2.4 Cerebellar hypoplasia
- 2.5 Heterotopia
- 2.6 Holoprosencephaly
- 2.7 Hydrocephalus
- 2.8 Leukodystrophy
- 2.9 Lissencephaly

#### 3. Movement abnormality

- 3.1 Ataxia
- 3.2 Chorea
- 3.3 Dystonia
- 3.4 Parkinsonism

#### 4. Neuromuscular abnormality

- 4.1 Muscular hypotonia
- 4.2 Muscular hypertonia
- 4.3 Hyperreflexia
- 4.4 Spasticity

#### 5. Seizures

- 5.1 Febrile seizures
- 5.2 Focal seizures
- 5.3 Generalized seizures

#### 6. Others

- 6.1 Craniosynostosis
- 6.2 Dementia
- 6.3 Encephalopathy
- 6.4 Headache / Migraine
- 6.5 Macrocephaly
- 6.6 Microcephaly
- 6.7 Neuropathy
- 6.8 Stroke

### METABOLISM

- 1. Creatine kinase
- 2. Decreased plasma carnitine
- 3. Hyperalaninemia
- 4. Hypoglycemia
- 5. Increased CSF lactate
- 6. Increased serum pyruvate
- 7. Ketosis
- 8. Lactic acidosis
- 9. Organic aciduria

### EYE

- 1. Blepharospasm
- 2. Cataract
- 3. Coloboma
- 4. Glaucoma
- 5. Microphthalmos
- 6. Nystagmus
- 7. Ophthalmoplegia
- 8. Optic atrophy
- 9. Ptosis
- 10. Retinitis pigmentosa
- 11. Retinoblastoma
- 12. Strabismus
- 13. Visual impairment

### MOUTH, THROAT AND EAR

- 1. Abnormality of dental color
- 2. Cleft lip / palate
- 3. Conductive hearing impairment
- 4. External ear malformation
- 5. Hypodontia
- 6. Sensorineural hearing impairment

### SKIN, INTEGUMENT AND SKELETAL

#### 1. Skeletal

- 1.1 Abnormal limb morphology
- 1.2 Abnormal vertebral column
- 1.3 Joint hypermobility
- 1.4 Multiple joint contractures
- 1.5 Polydactyly
- 1.6 Scoliosis
- 1.7 Syndactyly
- 1.8 Talipes equinovarus

### 2. Skin and integument

- 2.1 Abnormal skin pigmentation
- 2.2 Abnormal hair
- 2.3 Abnormal nail
- 2.4 Hyperextensible skin
- 2.5 Ichthyosis

### CARDIOVASCULAR

- 1. Angioedema
- 2. Aortic dilatation
- 3. Arrhythmia
- 4. Coarctation of aorta
- 5. Defect of atrial septum
- 6. Defect of ventricular septum
- 7. Dilated cardiomyopathy
- 8. Hypertrophic cardiomyopathy
- 9. Lymphedema
- 10. Malf. of heart and great vessels
- 11. Myocardial infarction
- 12. Tetralogy of Fallot

### GASTROINTESTINAL, GENITOURINARY, ENDOCRINE

#### 1. Gastrointestinal

- 1.1 Aganglionic megacolon
- 1.2 Constipation
- 1.3 Diarrhea
- 1.4 High hepatic transaminases
- 1.5 Gastroschisis
- 1.6 Hepatic failure
- 1.7 Hepatomegaly
- 1.8 Obesity
- 1.9 Pyloric stenosis
- 1.10 Vomiting

#### 2. Genitourinary

- 2.1 Hydronephrosis
- 2.2 Renal agenesis /hypoplasia
- 2.3 Renal cyst
- 2.4 Renal tubular dysfunction

### 3. Endocrine

- 3.1 Diabetes mellitus
- 3.2 Hypothyroidism
- 3.3 Hyperparathyroidism
- 3.4 Hypoparathyroidism
- 3.5 Hyperthyroidism

### REPRODUCTION

- 1. Abnormal external genitalia
- 2. Abnormal internal genitalia
- 3. Hypogonadism
- 4. Hypospadias
- 5. Infertility

### ONCOLOGY

- 1. Adenomatous polyposis
- 2. Breast carcinoma
- 3. Colorectal carcinoma
- 4. Leukemia
- 5. Myelofibrosis
- 6. Neoplasm of the lung
- 7. Neoplasm of the skin
- 8. Paraganglioma
- 9. Pheochromocytoma

### HEMATOLOGY AND IMMUNOLOGY

- 1. Abnormality of coagulation
- 2. Anemia
- 3. Immunodeficiency
- 4. Neutropenia
- 5. Pancytopenia
- 6. Abnormal hemoglobin
- 7. Splenomegaly
- 8. Thrombocytopenia

### PRENATAL AND DEVELOPMENT

- 1. Failure to thrive
- 2. Hemihypertrophy
- 3. Hydrops fetalis
- 4. IUGR
- 5. Oligohydramnios
- 6. Overgrowth
- 7. Polyhydramnios
- 8. Premature birth
- 9. Disproportionate short stature
- 10. Proportionate short stature
- 11. Tall stature

**OTHER** (INCLUDING DYSMORPHIC FACIAL FEATURES AND OTHER DESCRIPTORS):

Associated Condition(s)	Test Type	Test Name	Test Code	Sample Type
<b>AMINO ACID, ORGANIC ACID, FATTY ACID OXIDATION DISORDERS</b>				
Multiple	Biochemical Assay	Acylcarnitine Profile	B0210	DBS, WB, gDNA
Multiple	Biochemical Assay	Amino Acid Profile	B2020	DBS, WB, gDNA
2,4 Dienoyl-CoA Reductase Deficiency (DE RED)	Full Gene Analysis	<i>NADK2</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
2-methylbutyryl Glycinuria	Full Gene Analysis	<i>ACADSB</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
3-methylcrotonyl-CoA Carboxylase Deficiency (3-MCC Deficiency)	Targeted Variant Testing	3-MCC Deficiency Mutation Panel	D0410	DBS
3-methylglutaconic Aciduria, Type I	Full Gene Analysis	<i>AUH</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Argininemia	Full Gene Analysis	<i>ARG1</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Argininosuccinic Aciduria	Full Gene Analysis	<i>ASL</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Beta-ketothiolase Deficiency	Full Gene Analysis	<i>ACAT1</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Carnitine Palmitoyltransferase I Deficiency	Full Gene Analysis	<i>CPT1A</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Carnitine Palmitoyltransferase II Deficiency	Full Gene Analysis	<i>CPT2</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Carnitine Uptake Defect (CUD)	Full Gene Analysis	<i>SLC22A5</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Carnitine-acylcarnitine Translocase (CACT) Deficiency	Full Gene Analysis	<i>SLC25A20</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Citrullinemia Type I	Full Gene Analysis	<i>ASS1</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Citrullinemia Type II	Full Gene Analysis	<i>SLC25A13</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Cobalamin C Deficiency	Full Gene Analysis	<i>MMACHC</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Cobalamin D Deficiency	Full Gene Analysis	<i>MMADHC</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Glutaric Acidemia Type I	Targeted Variant Testing	Glutaric Acidemia Type I Mutation Panel	D0406	DBS
Glutaricaciduria, Type I	Full Gene Analysis	<i>GCDH</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
HMG-CoA Lyase Deficiency	Full Gene Analysis	<i>HMGCL</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Homocystinuria	Full Gene Analysis	<i>CBS</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Hypermethioninemia	Full Gene Analysis	<i>ADK</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Isobutyryl-CoA Dehydrogenase Deficiency	Full Gene Analysis	<i>ACAD8</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Isovaleric Acidemia	Targeted Variant Testing	Isovaleric Acidemia Mutation Panel	D0409	DBS
Isovaleric Acidemia	Full Gene Analysis	<i>IVD</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Long-chain 3-hydroxyacyl-CoA Dehydrogenase Deficiency (LCHADD)	Targeted Variant Testing	<i>LCHADD</i> Mutation Panel	D0407	DBS
Maple Syrup Urine Disease	Targeted Variant Testing	Maple Syrup Urine Disease Mutation Panel	D0401	DBS
Medium-chain Acyl-CoA Dehydrogenase Deficiency (MCADD)	Targeted Variant Testing	<i>MCADD</i> Mutation Panel	D0400	DBS
Medium-chain Acyl-CoA Dehydrogenase Deficiency (MCADD)	Full Gene Analysis	<i>ACADM</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Methylmalonic Acidemia	Targeted Variant Testing	Methylmalonic Acidemia Mutation Panel	D0411	DBS
Methylmalonic Acidemia	Full Gene Analysis	<i>MUT</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Multiple Carboxylase Deficiency	Full Gene Analysis	<i>HLCS</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Multiple Sulfatase Deficiency	Full Gene Analysis	<i>SUMF1</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Phenylketonuria (PKU)	Biochemical Assay	PKU Monitoring - Phenylalanine	B0018	DBS, WB
Phenylketonuria (PKU)	Full Gene Analysis	<i>PAH</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Propionic Acidemia	Targeted Variant Testing	Propionic Acidemia Mutation Panel	D0412	DBS
Short Chain 3-hydroxyacyl-CoA Dehydrogenase Deficiency (M/SCHADD)	Full Gene Analysis	<i>HADH</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Short-chain Acyl-CoA Dehydrogenase Deficiency (SCADD)	Full Gene Analysis	<i>ACADS</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA

Associated Condition(s)	Test Type	Test Name	Test Code	Sample Type
Tyrosinemia	Biochemical Assay	Tyrosinemia Monitoring - Succinylacetone and Tyrosine	B0022	DBS, WB
Tyrosinemia Type I	Full Gene Analysis	<i>FAH</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Tyrosinemia Type I	Biochemical	Succinylacetone (SUAC)	B0021	DBS, WB, gDNA
Tyrosinemia Type II	Full Gene Analysis	<i>TAT</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Tyrosinemia Type III	Full Gene Analysis	<i>HPD</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Very Long-chain Acyl-CoA Dehydrogenase Deficiency (VLCADD)	Full Gene Analysis	<i>ACADVL</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
<b>BIOTINIDASE DEFICIENCY</b>				
Biotinidase Deficiency	Biochemical Assay	Biotinidase Deficiency (Complete/Partial) - Biotinidase Deficiency Enzyme Analysis	B0001	DBS
Biotinidase Deficiency	Targeted Variant Testing	Biotinidase Deficiency Mutation Panel	D0402	DBS
Biotinidase Deficiency	Full Gene Analysis	<i>BTD</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
<b>CYSTIC FIBROSIS</b>				
Cystic Fibrosis	Biochemical Assay	IRT Analysis (Not valid after 90 days of age)	B0005	DBS
Cystic Fibrosis	Targeted Variant Testing	Cystic Fibrosis Mutation Panel	D3100	DBS
Cystic Fibrosis	Full Gene Analysis	<i>CFTR</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
<b>DUCHENNE MUSCULAR DYSTROPHY</b>				
Duchenne Muscular Dystrophy (DMD)	Biochemical Assay	Duchenne Muscular Dystrophy Creatine Kinase Activity	B0006	DBS
Duchenne Muscular Dystrophy (DMD)	Full Gene Analysis	<i>DMD</i> Gene Sequencing and Del/Dup Testing	D4045	DBS, WB, SV, gDNA
Duchenne Muscular Dystrophy (DMD)	Deletion/Duplication Analysis	<i>DMD</i> Del/Dup Testing	D5125	DBS, WB, SV, gDNA
<b>FRIEDREICH'S ATAXIA</b>				
Friedreich's Ataxia	Tandem Repeat Analysis	<i>FXN</i> Repeat Analysis	D5133	DBS, WB, gDNA
<b>GALACTOSEMIA</b>				
Galactosemia	Biochemical Assay	Galactosemia Monitoring - Galactose-1-phosphate uridylyltransferase Enzyme Analysis and Total Galactose	B0009	DBS
Galactosemia	Targeted Variant Testing	Galactosemia Mutation Panel	D0405	DBS
Galactosemia	Full Gene Analysis	<i>GALT</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Galactosemipimerase Deficiency	Full Gene Analysis	<i>GALE</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Galactokinase Deficiency	Full Gene Analysis	<i>GALK</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
<b>GLUCOSE-6-PHOSPHATE DEHYDROGENASE DEFICIENCY</b>				
Glucose-6-phosphate Dehydrogenase Deficiency	Biochemical Assay	Glucose-6-phosphate Dehydrogenase Deficiency (screening only)	B0011	DBS
Glucose-6-phosphate Dehydrogenase Deficiency	Targeted Variant Testing	Glucose-6-phosphate Dehydrogenase Deficiency Mutation Panel	D0404	DBS
Glucose-6-phosphate Dehydrogenase Deficiency	Full Gene Analysis	<i>G6PD</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
<b>LYSOSOMAL STORAGE DISORDERS - TESTING OPTIONS</b>				
Lysosomal Storage Disorders	Biochemical Assay	Lysosomal Storage Disease Enzyme Panel	B2040	DBS, WB
Lysosomal Storage Disorders	Full Gene Analysis	Lysosomal Storage Disorder Gene Sequencing Panel (12 Genes)	D3001	DBS, WB, SV, gDNA
Fabry Disease	Biochemical Assay	Alpha-Galactosidase A Enzyme Analysis	B0007	DBS, WB
Fabry Disease	Biochemical Assay	Globotriaosylsphingosine (lyso-Gb3) Monitoring	B0029	DBS, WB
Fabry Disease	Full Gene Analysis	<i>GLA</i> Gene Sequencing	D5033	DBS, WB, SV, gDNA
Gaucher Disease	Biochemical Assay	Glucocerebrosidase (Glucosylceramidase) Enzyme Analysis	B0010	DBS, WB
Gaucher Disease	Biochemical Assay	Glucosylsphingosine (lyso-Gb1) Monitoring	B0030	DBS, WB
Gaucher Disease	Full Gene Analysis	<i>GBA</i> Gene Sequencing	D5032	DBS, WB, SV, gDNA
Krabbe Disease	Biochemical Assay	Galactocerebrosidase Enzyme Analysis	B0012	DBS, WB

Associated Condition(s)	Test Type	Test Name	Test Code	Sample Type
Krabbe Disease	Biochemical Assay	Psychosine Biochemical Assay	B0028	DBS, WB
Krabbe Disease	Full Gene Analysis	<i>GALC</i> Gene Sequencing	D5031	DBS, WB, SV, gDNA
MPS I (Hurler Syndrome)	Biochemical Assay	Alpha-L-Iduronidase Enzyme Analysis	B0013	DBS, WB
MPS I (Hurler Syndrome)	Full Gene Analysis	<i>IDUA</i> Gene Sequencing	D5041	DBS, WB, SV, gDNA
MPS II (Hunter Syndrome)	Biochemical Assay	Iduronate 2-Sulfatase Enzyme Analysis	B0014	DBS, WB
MPS II (Hunter Syndrome)	Full Gene Analysis	<i>IDS</i> Gene Sequencing	D5042	DBS, WB, SV, gDNA
MPS IVA (Morquio A Syndrome)	Biochemical Assay	Galactosamine-6-Sulfatase Enzyme Analysis	B0015	DBS, WB
MPS IVA (Morquio A Syndrome)	Full Gene Analysis	<i>GALNS</i> Gene Sequencing	D5028	DBS, WB, SV, gDNA
MPS IVB (GM1 Gangliosidosis)	Biochemical Assay	$\beta$ -galactosidase Enzyme Analysis	B0025	DBS, WB
MPS IVB (GM1 Gangliosidosis)	Full Gene Analysis	<i>GLB1</i> Gene Sequencing	D5034	DBS, WB, SV, gDNA
MPS VI (Maroteaux-Lamy Syndrome)	Biochemical Assay	Arylsulfatase B Enzyme Analysis	B0016	DBS, WB
MPS VI (Maroteaux-Lamy Syndrome)	Full Gene Analysis	<i>ARSB</i> Gene Sequencing	D5009	DBS, WB, SV, gDNA
MPS VII (Sly Syndrome)	Biochemical Assay	$\beta$ -glucuronidase Enzyme Analysis	B0026	DBS, WB
Mucopolysaccharidosis VII	Full Gene Analysis	<i>GUSB</i> Gene Sequencing	D5035	DBS, WB, SV, gDNA
Multiple Sulfatase Deficiency	Full Gene Analysis	<i>SUMF1</i> Gene Sequencing	D5058	DBS, WB, SV, gDNA
Niemann Pick Disease Types A and B	Biochemical Assay	ACID Sphingomyelinase Enzyme Analysis	B0017	DBS, WB
Niemann Pick Disease Types A and B	Full Gene Analysis	<i>SMPD1</i> Gene Sequencing	D5057	DBS, WB, SV, gDNA
Pompe Disease	Biochemical Assay	ACID Alpha-Glucosidase Enzyme Analysis	B0019	DBS, WB
Pompe Disease	Full Gene Analysis	<i>GAA</i> Gene Sequencing	D5025	DBS, WB, SV, gDNA
Neuronal Ceroid Lipofuscinosis 2 (CLN2)	Biochemical Assay	Tripeptidyl peptidase 1 Enzyme Analysis	B0027	DBS, WB
Neuronal Ceroid Lipofuscinosis 2 (CLN2)	Full Gene Analysis	<i>TPP1</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
<b>SEVERE COMBINED IMMUNODEFICIENCY</b>				
Severe Combined Immunodeficiency (SCID)	Molecular DNA Screen	TREC Assay	D0416	DBS
Severe Combined Immunodeficiency (SCID)	Full Gene Analysis	<i>ADA</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Severe Combined Immunodeficiency (SCID)	Full Gene Analysis	<i>AK2</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Severe Combined Immunodeficiency (SCID)	Full Gene Analysis	<i>ATM</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Severe Combined Immunodeficiency (SCID)	Full Gene Analysis	<i>CD3D</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Severe Combined Immunodeficiency (SCID)	Full Gene Analysis	<i>CD3E</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Severe Combined Immunodeficiency (SCID)	Full Gene Analysis	<i>CD3Z</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Severe Combined Immunodeficiency (SCID)	Full Gene Analysis	<i>CORO1A</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Severe Combined Immunodeficiency (SCID)	Full Gene Analysis	<i>DCLRE1C</i> (Artemis) Gene Sequencing	D3100	DBS, WB, SV, gDNA
Severe Combined Immunodeficiency (SCID)	Full Gene Analysis	<i>DOCK8</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Severe Combined Immunodeficiency (SCID)	Full Gene Analysis	<i>FOXN1</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Severe Combined Immunodeficiency (SCID)	Full Gene Analysis	<i>IL2RG</i> SGene Sequencing	D3100	DBS, WB, SV, gDNA
Severe Combined Immunodeficiency (SCID)	Full Gene Analysis	<i>IL7R</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Severe Combined Immunodeficiency (SCID)	Full Gene Analysis	<i>JAK3</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Severe Combined Immunodeficiency (SCID)	Full Gene Analysis	<i>LIG4</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Severe Combined Immunodeficiency (SCID)	Full Gene Analysis	<i>NHEJ1</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Severe Combined Immunodeficiency (SCID)	Full Gene Analysis	<i>ORAI1</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA

Associated Condition(s)	Test Type	Test Name	Test Code	Sample Type
Severe Combined Immunodeficiency (SCID)	Full Gene Analysis	<i>PNP</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Severe Combined Immunodeficiency (SCID)	Full Gene Analysis	<i>PRKDC</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Severe Combined Immunodeficiency (SCID)	Full Gene Analysis	<i>PTPRC</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Severe Combined Immunodeficiency (SCID)	Full Gene Analysis	<i>RAC2</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Severe Combined Immunodeficiency (SCID)	Full Gene Analysis	<i>RAG1</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Severe Combined Immunodeficiency (SCID)	Full Gene Analysis	<i>RAG2</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Severe Combined Immunodeficiency (SCID)	Full Gene Analysis	<i>RMRP</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Severe Combined Immunodeficiency (SCID)	Full Gene Analysis	<i>STIM1</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Severe Combined Immunodeficiency (SCID)	Full Gene Analysis	<i>TBX1</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Severe Combined Immunodeficiency (SCID)	Full Gene Analysis	<i>ZAP70</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
<b>SICKLE CELL AND OTHER HEMOGLOBINOPATHIES</b>				
Sickle Cell and Other Hemoglobinopathies	Biochemical Assay	Isoelectric Focusing GEL Electrophoresis of Hemoglobiins	B0020	DBS
Sickle Cell and Other Hemoglobinopathies	Targeted Variant Testing	Sickle Cell and Other Hemoglobinopathies Mutation Panel	D0408	DBS
<b>SPINAL MUSCULAR ATROPHY (SMA)</b>				
Spinal Muscular Atrophy (SMA)	Deletion/Duplication Analysis	SMA Diagnostic Test	D5134	DBS, WB, gDNA
Spinal Muscular Atrophy (SMA)	Deletion/Duplication Analysis	SMA Carrier Screen	D5135	DBS, WB, gDNA
Spinal Muscular Atrophy (SMA)	Deletion/Duplication Analysis	<i>SMN2</i> Copy Number Test	D5136	DBS, WB, SV, gDNA
<b>OTHER</b>				
Congenital Adrenal Hyperplasia (CAH)	Biochemical Assay	Congenital adrenal hyperplasia - 17A Hydroxyprogesterone (17 OHP)	B0002	DBS
Congenital Adrenal Hyperplasia (CAH)	Full Gene Analysis	<i>CYP21A2</i> Gene Sequencing and Del/Dup Testing (by MLPA)	D5019	DBS, WB, SV, gDNA
Congenital Hypothyroidism	Biochemical Assay	Thyroid-Stimulating Hormone (TSH)	B0003	DBS
Congenital Hypothyroidism	Biochemical Assay	Thyroxine (T4)	B0004	DBS
Fragile X	Triplet Repeat Testing	<i>FMR1</i> Triplet Repeat (CGG) Testing	D4042	DBS, WB, SV, gDNA
X-linked Adrenoleukodystrophy	Biochemical Assay	X-Linked Adrenoleukodystrophy - C26:0 Lysophosphatidylcholine	B0023	DBS, WB
X-linked Adrenoleukodystrophy	Full Gene Analysis	<i>ABCD1</i> Gene Sequencing	D3100	DBS, WB, SV, gDNA
Multiple	Biochemical Assay	Post Mortem - Includes: 17-Hydroxyprogesterone, Acylcarnitines, Galactose, and <i>TSH</i>	B0024	DBS

DBS = Dried Blood Spots, WB = Whole Blood, SV = Saliva Swab, gDNA = Genomic DNA

The purpose of this Informed Consent Form (ICF) is to provide you with a description of the Test ordered, known risks and benefits of the Test, anonymization of personal health information ("PHI"), sample and data retention, research opportunities, and the reporting of secondary findings, if applicable.

## TEST INFORMATION

Your healthcare provider ("HCP") has recommended that you, or your dependent, receive clinical testing ("Test") indicated on the submitted Test Requisition Form ("Requisition"). For more information on the reasons your HCP has ordered the Test, and the disorders your HCP is testing for, please consult with your HCP. Providing a Sample and undergoing the Test is voluntary.

*Enzyme/Biomarker Test:* This type of test measures the presence or absence of enzymes/biomarkers and/or their level of activity in an individual. Only the enzymes/biomarkers identified on the requisition will be tested. Results from this type of Test may indicate the presence of a specific condition or conditions, and follow-up confirmatory testing may be recommended.

*Genetic/Genomic Test:* This type of Test analyzes one or more segments of your DNA depending on the assay requested. This Test is used to identify what, if any, DNA variant(s) you or your dependent possesses that is causing the specific disease, condition or risk you are being tested for. Identifying the mutation may be useful for diagnostic and treatment purposes, and allows at-risk family members to be tested. In some cases, we may not be able to determine with certainty which gene is actually causing a disease.

## TEST METHOD

If you consent to the Test, your HCP will take a sample of your and/or your dependent's blood, saliva, body fluid, tissue or other sample type. The Sample will be sent to Revvity's laboratories in the United States for the Test; the majority of testing will be performed at our laboratory headquarters in Pittsburgh, PA.

Under some circumstances, including inadequate or poor quality sample, an additional Sample may be required for Tests to be performed.

## TEST RESULTS

Your treating HCP has sole responsibility for all decisions concerning the possible management of your diagnosis and disease; Revvity will not provide a diagnosis. Revvity will report Test results via secure email, a secure internet portal, or fax. Your HCP is responsible for communicating with you regarding the results of the Test and may refer you or your dependent to a specialist for further clinical evaluation and confirmation of diagnosis, if applicable. Possible results for Genetic/Genomic Tests include:

1. *Positive:* A positive genetic test result may indicate that you are a carrier of, predisposed to, or have the specific disease or condition being tested for.
2. *Negative:* A negative result indicates that no disease-causing variant was identified in the Test performed. No Test can rule out all genetic diseases or conditions. A negative result does not guarantee that you are free from genetic disorders or other medical conditions.
3. *Inconclusive/Variant of Uncertain Significance:* A variant of uncertain significance (VOUS) result indicates that a DNA change was detected, but it is currently unknown if the variant is associated with a genetic disorder. A VOUS is not the same as a positive result and does not clarify whether there is an increased risk to develop a genetic disorder. The variant could be a benign change or it could be indicative of disease/disease-causing.
4. *Unexpected Results:* This Test may reveal an important genetic change that is not directly related to the reason for ordering this test. This information would be disclosed to your HCP if it potentially impacts medical care, and you have consented to receive this type of result.

## TEST REPORT

Variants are described as pathogenic, likely pathogenic, or variant of uncertain significance in genes interpreted to be responsible for, or potentially contributing to, a disease or condition. For testing performed on prenatal samples, or for screening of apparently healthy individuals, only variants classified as pathogenic or likely pathogenic will be reported.

When Whole Exome Sequencing (WES) or Whole Genome Sequencing (WGS) tests are ordered by your HCP, you may have the option to receive some findings not directly related to the reason for ordering the Test called "Secondary Findings". When Secondary Findings are requested, only Pathogenic or Likely Pathogenic findings will be reported, where applicable. Please read the Secondary Findings sections on page 3 and/or 4 of this consent form for more information. Secondary findings are not available for all Tests.

## INFORMATION ABOUT PARENTAL AND FAMILIAL SAMPLES

In some circumstances, it may be helpful for additional family members to undergo testing in order to provide information that can aid in the interpretation of the Test results. These Tests could be part of a WES/WGS TRIO Test or as stand-alone targeted testing. If the HCP recommends testing for additional family members, family members may have the option to receive information about secondary findings either as a part of the proband report or as a standalone parental report. A full analysis of the parental samples for secondary findings will only be completed if standalone reports are ordered (for an additional charge). In conjunction with proband testing, any variants reported in the proband will include inheritance information from family members included in the study.

## TEST LIMITATIONS

Due to current limitations in technology and incomplete knowledge of diseases and genes, some variants may not be detected by the Test ordered. There is a possibility that the Test result is uninterpretable or deemed of unknown significance. Further testing may be required when more information is gained. In rare circumstances, Test results may be suggestive of a condition different from that which was originally considered for the purpose of consenting to this Test. The Test may also find variants or genes that lead to conditions for which you currently do not have symptoms or may not be related to your current condition.

## TEST RISKS

Patients and family members may experience anxiety before, during, and/or after testing. Testing multiple family members may reveal that familial relationships are not biologically what they were assumed to be. For example, the Test may indicate non-paternity (the stated father of an individual is not the biological father) or consanguinity (the parents of an individual are closely related by blood). These biological relationships may need to be reported to the HCP who ordered the test.

Taking a blood or tissue sample from you and/or your child may lead to mild pain, bruising, swelling, redness, and a slight risk of infection. Light-headedness, fainting or nausea may occur if your HCP collects blood or tissue samples. These side-effects are typically brief and transient, but you should contact your HCP if you and/or your child require treatment. Under some circumstances an additional sample may be required for Tests to be performed.

A positive test result may limit your access to health insurance or life assurance coverage; for example, a life insurance company might ask you to provide genetic information indicating a disorder if this information is available to you. Please refer to information on the Genetic Information Nondiscrimination Act (GINA) and applicable local laws for more information.



## CONFIDENTIALITY

You have the right to confidential treatment of the Sample and your PHI. Your HCP will provide Revvity with Personal Health Information (“PHI”) such as your name, date of birth, gender and clinical symptoms to help track your sample and report results. To maintain confidentiality, the test results will only be released to the referring health care provider, to others designated by the referring health care provider as being involved in your care, to the ordering laboratory, to the patient/guardian, to other health care providers involved in your diagnosis and treatment, or as otherwise required by law or regulation. Unless required by law, Revvity will not disclose your PHI to any person or entity except with your written consent.

You and your HCP can control how your Sample and PHI are processed. You have the right to request access to your PHI, request corrections of any errors in recorded PHI, or where PHI may be missing or incomplete ask that it be completed. You also have the right to ask that your PHI be erased, subject to law or regulation. You can contact your HCP for such requests and your HCP will contact Revvity, or you can contact Revvity directly by visiting [www.revvity.com](http://www.revvity.com). If requests for access, correction, completion, or erasure cannot be fulfilled, you will be informed and provided with the reasons why your requests cannot be fulfilled.

## SAMPLE AND DATA RETENTION

Pursuant to laboratory best practices, your DNA sample will be retained by Revvity for a minimum of two years and then destroyed. Additionally, your PHI, the data from the Tests (including those performed before any withdrawal of consent) and the related reports will be retained by Revvity indefinitely, unless otherwise noted. In some instances, it may be beneficial to you for Revvity to retain your sample for a longer period of time in order to conduct additional testing, and Revvity will do so with appropriate documentation from you or your HCP.

Revvity is requesting consent to keep your and/or your child’s anonymized sample and data indefinitely for ongoing test development, scientific research, and/or other activities. Consent is optional, and the Test will be performed whether or not you provide consent to the following, as selected by your HCP on page 2 of the Test Requisition Form:

- Revvity will anonymize and retain your Sample indefinitely for internal quality control, test validation, assay development and improvement. By allowing Revvity to retain your Sample, you understand and agree that you give up any property rights you may have in the Sample and are donating it to Revvity Omics, Inc. If you withdraw your consent, no additional tests or anonymization will be carried out on your Sample; no results will be reported and your sample, reports and data that have not been anonymized will be destroyed.
- Revvity will anonymize your data and retain the anonymized data and related anonymized reports from your Tests indefinitely for internal statistical, quality analysis, research, scientific and technical development, and market research.

## RESEARCH OPTIONS

Revvity may collaborate with scientists, researchers and drug developers to advance knowledge of genetic diseases. If there are opportunities to participate in future research relevant to the disease in you and/or your child, Revvity may contact you or your HCP about the development of new testing, drug development, or other treatments. Revvity may also work with scientists or researchers from academic or commercial institutions who have received the necessary approvals to conduct a research study. In some instances, these scientists or researchers may like to contact you directly about your interest in participating in a specific research study. You may opt-out of research on page 2 of the Test Requisition Form.

## WITHDRAWAL OF CONSENT

I understand this consent is voluntary and is valid until I withdraw my consent. I understand I may withdraw my consent to sample and data retention, and to the Test at any time, that Revvity will not perform the Test unless consent has been obtained by the HCP. If I withdraw any consent, it will not affect actions taken before I withdrew my consent, including any anonymization of data or of my Sample. I understand that if I wish to withdraw my consent I should contact Revvity via email at: [genomics@revvity.com](mailto:genomics@revvity.com) or toll-free by telephone +1-866-354-2910 to request withdrawal.

## SECONDARY FINDINGS: APPLIES ONLY TO WES/WGS

- 1. ACMG Recommended Secondary Findings:** The American College of Medical Genetics and Genomics (ACMG), has recommended that secondary findings should be offered for a specific subset of highly penetrant and medically actionable genes associated with various inherited disorders for all individuals undergoing WGS or WES. Please refer to the latest version of the ACMG Recommendations for Reporting of Secondary Findings in Clinical Exome and Genome Sequencing for complete details of genes and conditions at [www.acmg.net](http://www.acmg.net). Medically-actionable conditions are those for which there is currently recommended treatment or preventative actions that can be taken to reduce the risk of developing the disease. An example would be hereditary cancer syndromes such as Lynch syndrome. Please note that only Pathogenic or Likely Pathogenic variants will be reported if this category of Secondary Findings is selected. We are unable to guarantee that the Test will find all medically-actionable conditions for which you have a pathogenic or likely pathogenic variant. You may have a pathogenic or likely pathogenic variant for a condition in which there was little or no coverage in the Test and therefore will not be detected. Additional testing for health purposes should be discussed with your doctor or genetic counselor.
- 2. Pharmacogenetic variants:** This category of Secondary Findings will include changes in the DNA that do not cause a disease but may be related to how your body processes certain medications, such as chemotherapy drugs, antipyretics, antidepressants, anticoagulants, and others. These variants may tell you how well medications will work or if you will have side effects if you do take the medications now or in the future.
- 3. Carrier status (ex. cystic fibrosis):** This category of Secondary Findings will include carrier findings for a select list of autosomal recessive conditions. The list of genes included in the carrier reporting is available upon request. A recessive condition is one in which two disease-causing variants in the same gene are typically required in order to show symptoms of the disease (one variant is inherited from each parent). Someone who has only one disease-causing variant is called a carrier. Please note that only Pathogenic or Likely Pathogenic variants will be reported if this category of Secondary Findings is selected. Further testing may be necessary to look for a second disease-causing variant in that gene. The Test is not designed to be a comprehensive carrier test. We are unable to guarantee that all conditions for which you are a carrier will be determined by the Test. You may be a carrier for a condition in which there was little or no coverage in the Testing and therefore will not be detected. Additional carrier testing for reproductive purposes should be discussed with your doctor or genetic counselor.
- 4. Diagnostic findings in all other disease-causing genes not related to your clinical features:** This category of Secondary Findings will include conditions that are medically-actionable but not included in the ACMG-recommended list, as well as conditions that are not medically-actionable (do not have recommended treatment or preventative measures). An example would be Alzheimer’s disease. Please note that only Pathogenic or Likely Pathogenic variants will be reported if this category of Secondary Findings is selected. Furthermore, we are unable to guarantee that the Test will find all disease-causing variants in all disease-causing genes. You may have a disease-causing variant for a condition in which there was little or no coverage in the Test and therefore will not be detected. Additional testing for health purposes should be discussed with your doctor or genetic counselor.