

Data Generation Requisition Form

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 Dried Blood Spots DNA, Source: Other:
 Collection Date: 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

SEND ADDITIONAL COPY OF RESULTS TO (If applicable)

Name Role with patient/Job title Clinic/Hospital/Institution Name

 Phone Number Fax Number Email Address

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				

Data Generation Requisition Form

STEP 4: TEST MENU

WHOLE EXOME SEQUENCING TESTING OPTIONS

D1500 Exome - Data Only (per sample)*

WHOLE GENOME SEQUENCING TESTING OPTIONS

D2500 Genome - Data Only (per sample)*

* FASTQ Data Files Only

DATA DELIVERY TYPE

Electronic Transfer

Hard Drive⁺

⁺ Additional Charges Apply

PERSON TO RECEIVE DATA

Name: _____

Street Address: _____

City/Town: _____ State: _____ Zip: _____

Phone: _____ Email: _____

STEP 5: DATA PERMISSIONS*

1. Check here if you would like to opt out of anonymized sample retention. Note, if not checked, this is interpreted as "consent given."
2. Check here if you would like to opt out of anonymized data retention. Note, if not checked, this is interpreted as "consent given."
3. Check here if you would like to opt out of your contact information being shared by Revvity to external researchers for direct communication regarding their studies. Note, if not checked, this is interpreted as "consent given."
4. Check here if sample(s) was or will be collected in the state of NV NY or OR. Additional consent must be completed, which can be found on the Revvity Omics website.

* For more information on data permissions options, please reference Page 4.

STEP 6: 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 _____