

STAT Prenatal and Fetal Testing 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: GESTATIONAL CARRIER'S INFORMATION (FETUS OF)

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

**Gestational Carrier's 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:**  Genomic DNA from fetal sample  Cultured Amniocytes  Cultured CVS  Direct Product of Conception (POC)  Fetal Tissue-Tissue Type: \_\_\_\_\_  Cord Blood  Other: \_\_\_\_\_ **contact lab before sending**  
 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				

## STAT Prenatal and Fetal Testing Requisition Form

### STEP 4: TEST MENU

#### STAT Whole Exome Sequencing Testing Options

- D1000E STAT Prenatal Whole Exome Sequencing, Proband ONLY
- D1330E STAT Prenatal Whole Exome Sequencing, DUO
- D1310E STAT Prenatal Whole Exome Sequencing, TRIO

#### STAT Whole Genome Sequencing Testing Options

- D0900E STAT Prenatal Prenatal CNGnome NGS Array
- D2010E STAT Prenatal Whole Genome Sequencing, Proband ONLY
- D2330E STAT Prenatal Whole Genome Sequencing, DUO
- D2310E STAT Prenatal Whole Genome Sequencing, TRIO

#### Additional Test Code Request (Non-Ongoing Pregnancies Only)

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

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

#### Targeted Testing Options

- D0600E STAT Prenatal Single-site testing

	MM / DD / YYYY
Proband Last Name, First Name	Proband DOB

Proband's Accession ID	Relationship to Proband

Positive Control Sample:  Already at Revvity  To be sent later  Not Available

Gene(s)	Coding Name (c.)	Protein Name (p.)

#### PREGNANCY STATUS

Ongoing pregnancy?  Yes  No

#### MATERNAL CELL CONTAMINATION (MCC) SAMPLE INFORMATION

For samples derived from CVS or amniotic fluid (AF), MCC is mandatory. For all other sample types, MCC is recommended but not required.

- D0990E MCC Test
- Submit MCC report performed in another lab

- D0999E STAT Prenatal Targeted CNV Analysis

	MM / DD / YYYY
Proband Last Name, First Name	Proband DOB

Proband's Accession ID	Relationship to Proband

Positive Control Sample:  Already at Revvity  To be sent later  Not Available

Cytoband/Gene	CN Event/Size/Exon	Additional CN Event/Size/Exon

**! Please include a copy of relative's report, if available.**

### STEP 5: FAMILIAL INFORMATION (REQUIRED WITH ALL TRIO/DUO AND MCC STUDIES)

#### FAMILY MEMBER 1

Last name, First name	Relationship to Patient

Date of Birth: MM / DD / YYYY    Symptomatic (clinically affected):  Yes  No    Sample  Included - Collection Date MM / DD / YYYY  To be sent later

Sample type:  Whole blood (preferred)  Saliva  Isolated DNA

#### FAMILY MEMBER 1

Last name, First name	Relationship to Patient

Date of Birth: MM / DD / YYYY    Symptomatic (clinically affected):  Yes  No    Sample  Included - Collection Date MM / DD / YYYY  To be sent later

Sample type:  Whole blood (preferred)  Saliva  Isolated DNA

**STAT Prenatal and Fetal Testing Requisition Form**

**STEP 6: PREGNANCY DETAILS AND FAMILY HISTORY**

**GESTATIONAL AGE AT SAMPLE COLLECTION**

Weeks: \_\_\_\_\_ Days: \_\_\_\_\_ EDD: \_\_\_\_\_ Based on:  LMP  Ultrasound

Gestation:  Singleton  Twins  Higher-Order Multiples

Was this pregnancy conceived by IVF?  Yes  No

Age of donor at egg retrieval: \_\_\_\_\_

Was an egg donor used to conceive this pregnancy?  Yes  No

Is the biological mother/egg donor the same as gestational carrier?  Yes  No

Fetal ultrasound abnormalities identified?

Yes  No

Fetal sex known?

Male  Female  Unknown  Ambiguous

Has gestational carrier had positive serum screen result?

Yes  No

Has biological mother/gamete donor had previous pregnancy with chromosome abnormality?

Yes  No

Has any previous genetic testing been performed on the fetus (including preimplantation testing on the embryo)?

Yes  No If Yes, what?

Does either biological parent or gamete donor have family history of:

• chromosome abnormality?  Yes  No If yes: \_\_\_\_\_

• genetic disease?  Yes  No If yes: \_\_\_\_\_

**STEP 7: 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 \_\_\_\_\_

**FETAL PHENOTYPES**

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

**PRIMARY INDICATION (required)**

- Fetal anomaly
- Cystic hygroma / Increased nuchal translucency size (mm): \_\_\_\_\_
- Abnormal NIPS result<sup>†</sup>
- Abnormal serum screen<sup>†</sup>
- Advanced maternal age<sup>†</sup>

<sup>†</sup> CNGnome® testing is recommended for this indication.

Whole genome and exome sequencing will not be performed without additional clinical information.

Previous test results should be sent with order.

**GASTROINTESTINAL**

- Absent stomach bubble
- Diaphragmatic hernia
- Echogenic bowel
- Gastroschisis
- Omphalocele

**GROWTH**

- Fetal demise
- Fetal pyelectasis/hydronephrosis
- Hydrops fetalis
- Intrauterine growth retardation (IUGR)
- Oligohydramnios
- Polyhydramnios
- Prematurity
- Sacrococcygeal teratoma
- Short long bones
- Small for gestational age (SGA)
- Small thorax

**NEUROLOGICAL**

- Abnormality of septum pellucidum
- Absent septum pellucidum
- Cavum septum pellucidum
- Choroid plexus cyst (CPC)
- Decreased fetal movement
- Encephalocele
- Myelomeningocele/Spina bifida

**CARDIAC**

- Congenital heart defect
- Intracardiac echogenic focus (IEF)
- Pericardial effusion

**OTHER**

- Absent nasal bone
- Fetal ascites
- Generalized edema
- Pleural effusion

Other: \_\_\_\_\_