



This requisition form can be used to submit a specimen for the Anemia Identified program, a no-charge US testing program sponsored by Agios and offered through Revvity. All patients suspected of having, or with a family history of, Hereditary Anemia (HA) can take part in the Anemia Identified Program.

The Anemia Identified program facilitates access to genetic testing to help in the diagnosis of HA or carrier status identification of HA. While Agios provides financial support for this program, tests and services are performed by an independent third party, Revvity. Healthcare providers must confirm that patients meet certain criteria to use the program. Agios receives de-identified patient data from this program, but at no time does Agios receive patient identifiable information. Agios receives contact information for healthcare providers who use this program. Genetic testing is available in the U.S. and Puerto Rico only. Healthcare providers who use this program have no obligation to recommend, purchase, order, prescribe, promote, administer, use or support any Agios product.

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: Saliva Swab Whole Blood Collection Date: Was this sample collected in the State of NY, NV or OR? Yes No

INDICATION FOR TESTING

Diagnosis in symptomatic patient Carrier testing Presymptomatic testing of at-risk family member Other: _____

STEP 2: ORDERING PROVIDER AND REPORTING PREFERENCE

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 Email Address

 Phone Number Fax Number

STEP 3: BILLING INFORMATION

INSTITUTIONAL BILLING

Agios B0374
 Institution/Organization Name Billing Account ID

Do not bill patient's insurance.

FOR INTERNAL USE ONLY				
Date Rec'd	TEMP	SPEC	COL	Rec'd #TUBES VOL



STEP 4: TEST MENU

- AGS001 Hereditary Anemia Panel
- AGS002 Pyruvate Kinase Enzyme Activity

Test Ordered	Acceptable Sample	Cold Pack Needed?
Hereditary Anemia Panel	1 Whole Blood (EDTA) sample or 1 saliva sample	NO
Pyruvate Kinase Enzyme Activity	1 Whole Blood (EDTA) sample	YES*
Hereditary Anemia Panel and Pyruvate Kinase Enzyme Activity together	2 Whole Blood (EDTA) samples	YES*

* Follow the instructions included in the whole blood collection pack to ensure proper usage of the cold pack.

STEP 5: PHYSICIAN CONFIRMATION OF INFORMED CONSENT AND MEDICAL NECESSITY

The undersigned person (or representative thereof) ensures he/she is a licensed medical professional authorized to order genetic testing and confirms that the patient has given appropriate informed consent for the testing ordered, including a discussion of the benefits and limitations. I confirm that testing is medically necessary and that test results may impact medical management for the patient. Furthermore, all information on this TRF is true to the best of my knowledge. My signature applies to the informed consent and/or attached letter of medical necessity.

I consent to Revvity sharing my name and contact information, and the date on which I executed this TRF, with Agios Pharmaceuticals and I understand and agree that I may be contacted by Agios Pharmaceuticals in connection with the genetic testing program for the purpose of receiving information on hereditary anemias, including pyruvate kinase deficiency.

Signature _____ Date _____

ADDITIONAL OPTIONAL PHENOTYPE / PATIENT HISTORY SECTION

Clinical History

CBC Data	Relevant Clinical Information
WBC: _____	<input type="radio"/> Asymptomatic <input type="radio"/> Symptomatic: _____
HGB: _____	<input type="radio"/> Acquired <input type="radio"/> Lifelong/familial <input type="radio"/> Perinatal/neonatal <input type="radio"/> Chronic <input type="radio"/> Episodic/sporadic
HCT: _____	Recent transfusion: <input type="radio"/> Yes <input type="radio"/> No Last transfusion date (mm-dd-yyyy): _____
RBC: _____	Family history: <input type="radio"/> Yes <input type="radio"/> No Disorder/relation to patient: _____
MCV: _____	Parental consanguinity: <input type="radio"/> Yes <input type="radio"/> No
MCH: _____	Blood smear shows: _____
CHC: _____	Bone marrow shows: _____
RDW: _____	
PLT: _____	
Rectics %: _____	
Abs Retic: _____	
Ferritin: _____	

Indication for Testing (See Metabolic Hematology Profile Comparison Chart)

Suspect	Previous Results
<input type="radio"/> Hereditary spherocytosis	Previous protein/functional testing: <input type="radio"/> Yes _____
<input type="radio"/> Hereditary elliptocytosis	Hb electrophoresis: _____
<input type="radio"/> Hereditary pyropoikilocytosis	
<input type="radio"/> Hereditary stomatocytosis	<input type="radio"/> G6PD activity level: _____ Coombs: <input type="radio"/> Pos <input type="radio"/> Neg <input type="radio"/> Not Done
<input type="radio"/> Pyruvate Kinase Deficiency	<input type="radio"/> IK activity level: _____ Splenectomy: <input type="radio"/> Yes <input type="radio"/> No
<input type="radio"/> Southeast Asian ovalocytosis	<input type="radio"/> Other enzyme level(s): _____
<input type="radio"/> Congenital dyserythropoietic anemia	<input type="radio"/> Osmotic fragility: <input type="radio"/> Normal <input type="radio"/> Increased <input type="radio"/> Decreased <input type="radio"/> Not performed
<input type="radio"/> Enzyme disorder: _____	<input type="radio"/> EMA binding/Band3: <input type="radio"/> Normal <input type="radio"/> Abnormal <input type="radio"/> Not performed
<input type="radio"/> Other: _____	<input type="radio"/> Ektacytometry: _____