

## Test Catalog for Pharma

Newborn Screening .....	2
Biochemical Genetic Testing .....	2
Molecular Genetic Testing .....	4
Genetic Testing .....	6
Molecular Genetic Testing .....	6
Biochemical Genetic Testing .....	10
How to order? .....	11

# Newborn Screening > Biochemical Genetic Testing

## StepOne® (Comprehensive with SCID)

### Fatty Acid Oxidation Disorders

Carnitine/Acylcarnitine Translocase Deficiency  
Carnitine Palmitoyl Transferase Deficiency Type I1  
3-Hydroxy Long Chain Acyl-CoA Dehydrogenase Deficiency  
2,4-Dienoyl-CoA Reductase Deficiency1  
Medium Chain Acyl-CoA Dehydrogenase Deficiency  
Multiple Acyl-CoA Dehydrogenase Deficiency  
Neonatal Carnitine Palmitoyl Transferase Deficiency Type II  
Short Chain Acyl-CoA Dehydrogenase Deficiency  
Short Chain Hydroxy Acyl-CoA Dehydrogenase Deficiency  
Trifunctional Protein Deficiency  
Very Long Chain Acyl-CoA Dehydrogenase Deficiency

### Organic Acid Disorders

3-Hydroxy-3-Methylglutaryl-CoA Lyase Deficiency  
Glutaric Acidemia Type I  
Isobutyryl-CoA Dehydrogenase Deficiency  
Isovaleric Acidemia  
2-Methylbutyryl-CoA Dehydrogenase Deficiency  
3-Methylcrotonyl-CoA Carboxylase Deficiency  
3-Methylglutaconyl-CoA Hydratase Deficiency  
Methylmalonic Acidemias  
Methylmalonyl-CoA Mutase Deficiency  
Some Adenosylcobalamin Synthesis Defects  
Maternal Vitamin B12 Deficiency  
Mitochondrial Acetoacetyl-CoA Thiolase Deficiency  
Propionic Acidemia  
Multiple CoA Carboxylase Deficiency  
Malonic Aciduria

### Amino Acid Disorders

Argininemia  
Argininosuccinic Aciduria  
5-Oxoprolinuria1  
Carbamoyl Phosphate Synthetase Deficiency1  
Citrullinemia  
Homocystinuria  
Hypermethioninemia  
Hyperammonemia, Hyperornithinemia, Homocitrullinuria Syndrome1  
Hyperornithinemia with Gyral Atrophy1  
Maple Syrup Urine Disease  
Phenylketonuria  
Classical/Hyperphenylalaninemia  
Biotin Cofactor Deficiencies  
Tyrosinemia  
Transient Neonatal Tyrosinemia  
Tyrosinemia Type I2  
Tyrosinemia Type II  
Tyrosinemia Type III

### Other Observations

Hyperalbuminemia  
Liver Disease  
Medium Chain Triglyceride Oil Administration  
Presence of EDTA Anticoagulants in blood specimen  
Treatment with Benzoate, Pyruvic Acid, or Valproic Acid  
Carnitine Uptake Deficiency

### Endocrine Disorder

Congenital Adrenal Hyperplasia  
Salt Wasting 21-Hydroxylase Deficiency  
Simple Virilizing 21-Hydroxylase Deficiency  
Congenital Hypothyroidism

### Other Disorder

Biotinidase Deficiency  
Complete Deficiency  
Partial Deficiency  
Glucose-6-Phosphate Dehydrogenase Deficiency  
Cystic Fibrosis (not valid after 3 months of age)\*  
Galactosemia  
Galactokinase Deficiency  
Galactose-1-Phosphate Uridyltransferase Deficiency  
Galactose-4-Epimerase Deficiency  
Severe Combined Immunodeficiency (SCID)

### Hemoglobin Disorder

Sickle Cell & other Hemoglobinopathies  
Hemoglobin S, S/C, S/Beta-Thalassemia, C, & E Diseases

## StepOne + SCID + LSDs

### StepOne (Comp + SCID)

(All disorders listed above)

### Lysosomal Storage Disorders (LSDs)

Fabry  
Gaucher  
Krabbe Disease  
Mucopolysaccharidosis Type I (MPS-I)  
Niemann-Pick A/B  
Pompe

## StepOne + LSDs + X-ALD (full panel)

### StepOne (Comp + SCID)

(All disorders listed above)

### Lysosomal Storage Disorders (LSDs)

Fabry  
Gaucher  
Krabbe Disease  
Mucopolysaccharidosis Type I (MPS-I)  
Niemann-Pick A/B  
Pompe

### X-Linked Adrenoleukodystrophy (X-ALD)

DMD

## Newborn Screening > Biochemical Genetic Testing

### LSDs Panel Only

---

#### Lysosomal Storage Disorders (LSDs)

Fabry  
Gaucher  
Krabbé Disease  
Mucopolysaccharidosis Type I (MPS-I)  
Niemann-Pick A/B  
Pompe

### StepOne® SCID only

---

### Specialty Testing

---

#### Post Mortem Screens

#### PKU Clinical Monitoring

#### Disclaimer

*The analyses conducted by PerkinElmer® Genetics produce results that can be used by qualified physicians in the diagnosis of disorders described herein. Evidence of these conditions will be detected in the vast majority of affected individuals; however, due to genetic variability, age of the patient at the time of specimen collection, quality of the specimen, health status of the patient, and other variables, such conditions may not be detected in all affected patients.*

*Pursuant to applicable federal and/or state laboratory requirements, PerkinElmer® Genetics has established and verified the accuracy and precision of its laboratory developed testing services. Tests are developed and performance characteristics are determined by PerkinElmer Genetics.*

# Newborn Screening > Molecular Genetic Testing

## Second Tier Reflex Testing Menu

PerkinElmer Genetics uses combinations of assays in a multi-tier approach that optimizes detection of abnormal results. Positive DNA identification for many disorders further speeds definitive diagnosis and implementation of critical therapies.

### Biochemical Second Tier Testing

Disorder	Testing Approach
Congenital Adrenal Hyperplasia	First Tier; 17-OH P Second Tier; Extracted 17-OH P on all elevated.
Congenital Hypothyroidism	First Tier; either T4 or TSH. Second Tier TSH with a primary T4.
Galactosemia	First Tier; Total Galactose plus quantitative Uridyltransferase. Second Tier; Fractionated Galactose.

### DNA Second Tier Testing

Disorder	Mutations Detected
Galactosemia	N314D (Duarte) Q188R, S135L, K285N, and L195P (Classical)
Hemoglobinopathies	Hb S (173A>T), Hb C (172G>A), Hb E (232G>A), Hb D (121G>C) and Hb O (121G>A) β Thalassemias -29A>G, -88C>T, and IVS1+6T>C
Cystic Fibrosis	This chart contains the 23 mutations recommended by the ACOG/ACMG:

ΔF508	1717-1G>A	W1282X	2307insA
ΔI507	R560T	N1303K	Y1092X
G542X	R553X	394delTT	M1101K
G85E	G551D	Y122X	S1255X
R117H	1898+1G>A	R347H	3876delA
621+1G>T	2184delA	V520F	3905insT
711+1G>T	2789+5G>A	A559T	5/7/9T
1078delT	3120+1G>A	S549N	F508C
R334W	R1162X	S549R	I507V
R347P	3659delC	1898+5G>T	I506V
A455E	3849+10kbC>T	2183AA>G	

# Newborn Screening > Molecular Genetic Testing

## Second Tier Reflex Testing Menu

---

Disorder	Mutations Detected
Biotinidase Deficiency	G98:d7i3, Q456H, R157H, R538C, D252G and D444H; D444H;A171T, D444H; F403V, D444H;R157H
MCAD	A985A>G, 199T>C
LCHAD	1528G>C
Glutaric Acidemia 1	A421V (Amish) R402W (Caucasian)
Propionic Acidemia	E168K (Spanish) 1218del14/ins12 (Caucasian) 1170insT
Methylmalonic Acidemia	N219Y (Caucasian) G717V (African American)
3-Methylcrotonyl-CoA Carboxylase Def.	518insT (Mennonite)
Maple Syrup Urine Disease	Y438N (previously known as Y393N)
Isovaleric Acidemia	A282V

### Disclaimer

The analyses conducted by PerkinElmer® Genetics produce results that can be used by qualified physicians in the diagnosis of disorders described herein. Evidence of these conditions will be detected in the vast majority of affected individuals; however, due to genetic variability, age of the patient at the time of specimen collection, quality of the specimen, health status of the patient, and other variables, such conditions may not be detected in all affected patients.

Pursuant to applicable federal and/or state laboratory requirements, PerkinElmer® Genetics has established and verified the accuracy and precision of its laboratory developed testing services. Tests are developed and performance characteristics are determined by PerkinElmer Genetics.

## Genetic Testing > Molecular Genetic Testing

### Clinical Whole Genome

---

Contact us on 1-866-463-6436 for details

### Clinical Exome

---

Contact us on 1-866-463-6436 for details

### NeoSeq (NGS)

---

Contact us on 1-866-463-6436 for details

### LSD NGS panel (single gene or full panel sequencing)

---

**Fabry *GLA* gene sequencing**

**Gaucher *GBA* gene sequencing**

**Pompe *GAA* gene sequencing**

**Krabbe *GALC* gene sequencing**

**Niemann-Pick A/B *SMPD1* gene sequencing**

**Hurler Syndrome (MPS-I) *IDUA* gene sequencing**

**Hunter syndrome (MPS-II) *IDS* gene sequencing**

**Morquio syndrome type A (MPS IVA) *GALNS* gene sequencing**

**Beta-1 Galactosidase (Morquio syndrome type B, MPS IVB) *GLB1* gene sequencing**

**Maroteaux-Lamy (MPS-VI) *ARSB* gene sequencing**

**Mucopolysaccharidosis VII *GUSB* gene sequencing**

#### Disclaimer

*The analyses conducted by PerkinElmer® Genetics produce results that can be used by qualified physicians in the diagnosis of disorders described herein. Evidence of these conditions will be detected in the vast majority of affected individuals; however, due to genetic variability, age of the patient at the time of specimen collection, quality of the specimen, health status of the patient, and other variables, such conditions may not be detected in all affected patients.*

*Pursuant to applicable federal and/or state laboratory requirements, PerkinElmer® Genetics has established and verified the accuracy and precision of its laboratory developed testing services. Tests are developed and performance characteristics are determined by PerkinElmer Genetics.*

## Genetic Testing > Molecular Genetic Testing

### Second Tier Reflex Testing Menu

PerkinElmer Genetics uses combinations of assays in a multi-tier approach that optimizes detection of abnormal results. Positive DNA identification for many disorders further speeds definitive diagnosis and implementation of critical therapies.

#### Biochemical Second Tier Testing

Disorder	Testing Approach
Congenital Adrenal Hyperplasia	First Tier; 17-OH P Second Tier; Extracted 17-OH P on all elevated.
Congenital Hypothyroidism	First Tier; either T4 or TSH. Second Tier TSH with a primary T4.
Galactosemia	First Tier; Total Galactose plus quantitative Uridyltransferase. Second Tier; Fractionated Galactose.

#### DNA Second Tier Testing

Disorder	Mutations Detected
Galactosemia	N314D (Duarte) Q188R, S135L, K285N, and L195P (Classical)
Hemoglobinopathies	Hb S (173A>T), Hb C (172G>A), Hb E (232G>A), Hb D (121G>C) and Hb O (121G>A) β Thalassemias -29A>G, -88C>T, and IVS1+6T>C
Cystic Fibrosis	This chart contains the 23 mutations recommended by the ACOG/ACMG:

ΔF508	1717-1G>A	W1282X	2307insA
ΔI507	R560T	N1303K	Y1092X
G542X	R553X	394delTT	M1101K
G85E	G551D	Y122X	S1255X
R117H	1898+1G>A	R347H	3876delA
621+1G>T	2184delA	V520F	3905insT
711+1G>T	2789+5G>A	A559T	5/7/9T
1078delT	3120+1G>A	S549N	F508C
R334W	R1162X	S549R	I507V
R347P	3659delC	1898+5G>T	I506V
A455E	3849+10kbC>T	2183AA>G	

## Genetic Testing > Molecular Genetic Testing

### Second Tier Reflex Testing Menu

---

Disorder	Mutations Detected
Biotinidase Deficiency	G98:d7i3, Q456H, R157H, R538C, D252G and D444H; D444H;A171T, D444H; F403V, D444H;R157H
MCAD	A985A>G, 199T>C
LCHAD	1528G>C
Glutaric Acidemia 1	A421V (Amish) R402W (Caucasian)
Propionic Acidemia	E168K (Spanish) 1218del14/ins12 (Caucasian) 1170insT
Methylmalonic Acidemia	N219Y (Caucasian) G717V (African American)
3-Methylcrotonyl-CoA Carboxylase Def.	518insT (Mennonite)
Maple Syrup Urine Disease	Y438N (previously known as Y393N)
Isovaleric Acidemia	A282V

#### Disclaimer

*The analyses conducted by PerkinElmer® Genetics produce results that can be used by qualified physicians in the diagnosis of disorders described herein. Evidence of these conditions will be detected in the vast majority of affected individuals; however, due to genetic variability, age of the patient at the time of specimen collection, quality of the specimen, health status of the patient, and other variables, such conditions may not be detected in all affected patients.*

*Pursuant to applicable federal and/or state laboratory requirements, PerkinElmer® Genetics has established and verified the accuracy and precision of its laboratory developed testing services. Tests are developed and performance characteristics are determined by PerkinElmer Genetics.*



## Genetic Testing > Molecular Genetic Testing

### Clinical Whole Genome

---

Contact us on 1-866-463-6436 for details

### Clinical Exome

---

Contact us on 1-866-463-6436 for details

### NeoSeq (NGS)

---

Contact us on 1-866-463-6436 for details

### LSD NGS panel (single gene or full panel sequencing)

---

**Fabry *GLA* gene sequencing**

**Gaucher *GBA* gene sequencing**

**Pompe *GAA* gene sequencing**

**Krabbe *GALC* gene sequencing**

**Niemann-Pick A/B *SMPD1* gene sequencing**

**Hurler Syndrome (MPS-I) *IDUA* gene sequencing**

**Hunter syndrome (MPS-II) *IDS* gene sequencing**

**Morquio syndrome type A (MPS IVA) *GALNS* gene sequencing**

**Beta-1 Galactosidase (Morquio syndrome type B, MPS IVB) *GLB1* gene sequencing**

**Maroteaux-Lamy (MPS-VI) *ARSB* gene sequencing**

**Mucopolysaccharidosis VII *GUSB* gene sequencing**

#### Disclaimer

*The analyses conducted by PerkinElmer® Genetics produce results that can be used by qualified physicians in the diagnosis of disorders described herein. Evidence of these conditions will be detected in the vast majority of affected individuals; however, due to genetic variability, age of the patient at the time of specimen collection, quality of the specimen, health status of the patient, and other variables, such conditions may not be detected in all affected patients.*

*Pursuant to applicable federal and/or state laboratory requirements, PerkinElmer® Genetics has established and verified the accuracy and precision of its laboratory developed testing services. Tests are developed and performance characteristics are determined by PerkinElmer Genetics.*

### StepOne® (Comp + SCID)

---

#### Fatty Acid Oxidation Disorders

Carnitine/Acylcarnitine Translocase Deficiency  
Carnitine Palmitoyl Transferase Deficiency Type I1  
3-Hydroxy Long Chain Acyl-CoA Dehydrogenase Deficiency  
2,4-Dienoyl-CoA Reductase Deficiency1  
Medium Chain Acyl-CoA Dehydrogenase Deficiency  
Multiple Acyl-CoA Dehydrogenase Deficiency  
Neonatal Carnitine Palmitoyl Transferase Deficiency Type II  
Short Chain Acyl-CoA Dehydrogenase Deficiency  
Short Chain Hydroxy Acyl-CoA Dehydrogenase Deficiency  
Trifunctional Protein Deficiency  
Very Long Chain Acyl-CoA Dehydrogenase Deficiency

#### Organic Acid Disorders

3-Hydroxy-3-Methylglutaryl-CoA Lyase Deficiency  
Glutaric Acidemia Type I  
Isobutyryl-CoA Dehydrogenase Deficiency  
Isovaleric Acidemia  
2-Methylbutyryl-CoA Dehydrogenase Deficiency  
3-Methylcrotonyl-CoA Carboxylase Deficiency  
3-Methylglutaconyl-CoA Hydratase Deficiency  
Methylmalonic Acidemias  
Methylmalonyl-CoA Mutase Deficiency  
Some Adenosylcobalamin Synthesis Defects  
Maternal Vitamin B12 Deficiency  
Mitochondrial Acetoacetyl-CoA Thiolase Deficiency  
Propionic Acidemia  
Multiple CoA Carboxylase Deficiency  
Malonic Aciduria

#### Amino Acid Disorders

Argininemia  
Argininosuccinic Aciduria  
5-Oxoprolinuria1  
Carbamoyl Phosphate Synthetase Deficiency1  
Citrullinemia  
Homocystinuria  
Hypermethioninemia  
Hyperammonemia, Hyperornithinemia, Homocitrullinuria Syndrome1  
Hyperornithinemia with Gyral Atrophy1  
Maple Syrup Urine Disease  
Phenylketonuria  
Classical/Hyperphenylalaninemia  
Biopterin Cofactor Deficiencies  
Tyrosinemia  
Transient Neonatal Tyrosinemia  
Tyrosinemia Type I2  
Tyrosinemia Type II  
Tyrosinemia Type III

#### Other Observations

Hyperalbuminemia  
Liver Disease  
Medium Chain Triglyceride Oil Administration  
Presence of EDTA Anticoagulants in blood specimen  
Treatment with Benzoate, Pyvalic Acid, or Valproic Acid  
Carnitine Uptake Deficiency

#### Endocrine Disorder

Congenital Adrenal Hyperplasia  
Salt Wasting 21-Hydroxylase Deficiency  
Simple Virilizing 21-Hydroxylase Deficiency  
Congenital Hypothyroidism

#### Other Disorder

Biotinidase Deficiency  
Complete Deficiency  
Partial Deficiency  
Glucose-6-Phosphate Dehydrogenase Deficiency  
Cystic Fibrosis (not valid after 3 months of age)\*  
Galactosemia  
Galactokinase Deficiency  
Galactose-1-Phosphate Uridyltransferase Deficiency  
Galactose-4-Epimerase Deficiency  
Severe Combined Immunodeficiency (SCID)

#### Hemoglobin Disorder

Sickle Cell & other Hemoglobinopathies  
Hemoglobin S, S/C, S/Beta-Thalassemia, C, & E Diseases

### X-ALD only

---

#### X-Linked Adrenoleukodystrophy (X-ALD)

#### Disclaimer

*The analyses conducted by PerkinElmer® Genetics produce results that can be used by qualified physicians in the diagnosis of disorders described herein. Evidence of these conditions will be detected in the vast majority of affected individuals; however, due to genetic variability, age of the patient at the time of specimen collection, quality of the specimen, health status of the patient, and other variables, such conditions may not be detected in all affected patients.*

*Pursuant to applicable federal and/or state laboratory requirements, PerkinElmer® Genetics has established and verified the accuracy and precision of its laboratory developed testing services. Tests are developed and performance characteristics are determined by PerkinElmer Genetics.*

## How to order?

### Ordering and Results

---

Please [contact](#) PerkinElmer representative.

#### Results

[Click here to visit Results center](#)

To obtain an ID to download results from our secure portal, please contact Client Services to obtain the appropriate form.

#### Disclaimer

*The analyses conducted by PerkinElmer® Genetics produce results that can be used by qualified physicians in the diagnosis of disorders described herein. Evidence of these conditions will be detected in the vast majority of affected individuals; however, due to genetic variability, age of the patient at the time of specimen collection, quality of the specimen, health status of the patient, and other variables, such conditions may not be detected in all affected patients.*

*Pursuant to applicable federal and/or state laboratory requirements, PerkinElmer® Genetics has established and verified the accuracy and precision of its laboratory developed testing services. Tests are developed and performance characteristics are determined by PerkinElmer Genetics.*