



Whole Exome and Whole Genome Sequencing Secondary Findings | ACMG Incidental Findings

Gene	Gene MIM	Disease/Phenotype	Inheritance	Typical age of onset
ACTA2	102620	Familial thoracic aortic aneurysm; Multisystemic smooth muscle dysfunction syndrome; Moyamoya disease	AD	Child / Adult
ACTC1	102540	ACTC1 - Related Disorders	AD	Child / Adult
ACVRL1	601284	Hereditary hemorrhagic telangiectasia	AD	Child / Adult
APC	611731	APC-Associated Polyposis Conditions	AD	Child / Adult
APOB	107730	Familial hypercholesterolemia; Hypobetalipoproteinemia	AD/AR	Child / Adult
ATP7B	606882	Wilson disease	AR	Child / Adult
BMPR1A	601299	Juvenile polyposis syndrome	AD	Child / Adult
BRCA1	113705	BRCA1 - related cancer susceptibility	AD	Adult
BRCA2	600185	BRCA2 - related cancer susceptibility; Fanconi Anemia	AD/AR	Adult
BTD	609019	Biotinidase deficiency	AR	Child / Adult
CACNA1S	114208	CACNA1S - Related Disorders	AD	Child / Adult
CASQ2	114251	Catecholaminergic polymorphic ventricular tachycardia	AR	Child / Adult
COL3A1	120180	Ehlers-Danlos syndrome, vascular type	AD	Child / Adult
DSC2	125645	Arrhythmogenic right ventricular cardiomyopathy	AD	Child / Adult
DSG2	125671	Arrhythmogenic right ventricular cardiomyopathy	AD	Child / Adult

Gene	Gene MIM	Disease/Phenotype	Inheritance	Typical age of onset
DSP	125647	DSP - Related Disorders	AD/AR	Child / Adult
ENG	131195	Hereditary hemorrhagic telangiectasia	AD	Child / Adult
FBN1	134797	FBN1 - Related Disorders	AD	Child / Adult
FLNC	102565	FLNC - Related Disorders	AD	Adult
GAA	606800	Pompe disease	AR	Child / Adult
GLA	300644	Fabry disease	XL	Child / Adult
HFE	613609	Hereditary hemochromatosis (c.845G>A; p.C282Y homozygotes only)	AR	Adult
HNF1A	142410	Maturity-Onset of Diabetes of the Young	AD	Child / Adult
KCNH2	152427	Long-QT syndrome type 2	AD	Child / Adult
KCNQ1	607542	KCNQ1 - Related Disorders	AD / AR	Child / Adult
LDLR	606945	Familial hypercholesterolemia	AD	Child / Adult
LMNA	150330	LMNA - Related Disorders	AD/AR	Child / Adult
MAX	154950	Hereditary paraganglioma- pheochromocytoma syndrome	AD	Child / Adult
MEN1	613733	Multiple endocrine neoplasia type 1	AD	Child / Adult
MLH1	120436	Lynch syndrome	AD	Adult
MSH2	609309	Lynch syndrome	AD	Adult
MSH6	600678	Lynch syndrome	AD	Adult
MUTYH	604933	MUTYH -associated polyposis	AR	Adult
MYBPC3	600958	Hypertrophic cardiomyopathy; Dilated cardiomyopathy, Left ventricular noncompaction	AD	Child / Adult
MYH11	160745	Familial thoracic aortic aneurysm	AD	Child / Adult
MYH7	160760	MYH7 - Related Disorders	AD	Child / Adult
MYL2	160781	Hypertrophic cardiomyopathy	AD	Child / Adult
MYL3	160790	Hypertrophic cardiomyopathy	AD	Child / Adult
NF2	607379	Neurofibromatosis type 2	AD	Child / Adult
OTC	300461	Ornithine transcarbamylase deficiency	XL	Child
PALB2	610355	Hereditary breast cancer; Fanconi Anemia	AD/AR	Adult
PCSK9	607786	Familial hypercholesterolemia	AD	Child / Adult
PKP2	602861	Arrhythmogenic right ventricular cardiomyopathy	AD	Child / Adult
PMS2	600259	Lynch syndrome; mismatch repair cancer syndrome	AD/AR	Adult

Gene	Gene MIM	Disease/Phenotype	Inheritance	Typical age of onset
PRKAG2	602743	Hypertrophic cardiomyopathy; Glycogen storage disease of heart; Wolff-Parkinson-White syndrome	AD	Child / Adult
PTEN	601728	PTEN hamartoma tumor syndrome	AD	Child / Adult
RB1	614041	Retinoblastoma	AD	Child / Adult
RET	164761	Multiple endocrine neoplasia type 2 A/B; Familial medullary thyroid carcinoma	AD	Child / Adult
RPE65	180069	RPE65 -related retinopathy	AR	Child / Adult
RYR1	180901	RYR1 - Related Disorders	AD/AR	Child / Adult
RYR2	180902	Catecholaminergic polymorphic ventricular tachycardia; Arrhythmogenic right ventricular dysplasia	AD	Child / Adult
SCN5A	600163	SCN5A - Related Disorders	AD	Child / Adult
SDHAF2	613019	Hereditary paraganglioma-pheochromocytoma syndrome	AD	Child / Adult
SDHB	185470	Hereditary paraganglioma-pheochromocytoma syndrome	AD	Child / Adult
SDHC	602413	Hereditary paraganglioma-pheochromocytoma syndrome	AD	Child / Adult
SDHD	602690	Hereditary paraganglioma-pheochromocytoma syndrome	AD	Child / Adult
SMAD3	603109	Loeys-Dietz syndrome	AD	Child / Adult
SMAD4	600993	SMAD4 - Related Disorders	AD	Child / Adult
STK11	602216	Peutz-Jeghers syndrome	AD	Child / Adult
TGFBR1	190181	Loeys-Dietz syndrome	AD	Child / Adult
TGFBR2	190182	Loeys-Dietz syndrome	AD	Child / Adult
TMEM127	613403	Hereditary paraganglioma-pheochromocytoma syndrome	AD	Child / Adult
TMEM43	612048	Arrhythmogenic right ventricular cardiomyopathy	AD	Child / Adult
TNNI3	191044	Hypertrophic cardiomyopathy	AD	Child / Adult
TNNT2	191045	TNNT2 - Related Disorders	AD	Child / Adult
TP53	191170	Li-Fraumeni syndrome	AD	Child / Adult
TPM1	191010	Hypertrophic cardiomyopathy; Dilated cardiomyopathy, Left ventricular noncompaction	AD	Child / Adult
TRDN	603283	Long QT syndrome; catecholaminergic polymorphic ventricular tachycardia	AR	Child / Adult

Gene	Gene MIM	Disease/Phenotype	Inheritance	Typical age of onset
TSC1	605284	Tuberous sclerosis complex	AD	Child
TSC2	191092	Tuberous sclerosis complex	AD	Child
TTN	188840	TTN - Related Disorders (DCM for ACMG)	AD/AR	Child / Adult
VHL	608537	Von Hippel-Lindau syndrome	AD	Child / Adult
WT1	607102	WT1 -related Wilms tumor	AD	Child
BAG3	603883	Dilated cardiomyopathy	AD	mainly adult, ped reported
DES	125660	Dilated cardiomyopathy	AD/AR	mainly adult, ped reported
RBM20	613171	Dilated cardiomyopathy	AD	Child / Adult
TNNC1	191040	Dilated cardiomyopathy	AD	mainly adult, ped reported
TTR	176300	Hereditary TTR amyloidosis	AD	Adult

How to order?

Please reach out to your nearest Revvity Omics' representatives to order for any queries, feedback, or information you may require. You may also contact us on:

- **US**

Email: genomics@revvity.com
Phone: 866-354-2910

- **UK**

Email: uk.genomics@revvity.com

- **Sweden**

Email: sweden.genomics@revvity.com

- **India**

Email: customercare.dxindia@revvity.com
Phone: 1800-266-6660 | 1800-266-1660

- **China**

Email: iclserviceg.china@revvity.com
Phone: 400-110-8718

Visit our website at www.revvityomics.com

Disclaimer: This testing service has not been cleared or approved by the U.S. Food and Drug Administration. Testing services may not be licensed in accordance with the laws in all countries. The availability of specific test offerings is dependent upon laboratory location. The content of this pamphlet is provided for informational purposes only, not as medical advice. It is not intended to substitute the consultation, diagnosis, and/or treatment provided by a qualified licensed physician or other medical professional.

