

Newborn sequencing research - reimagined.

Official gene panel

Introducing the fast-track to genetic discovery.

Revvity is powering easier and faster genetic interpretation for genetic and newborn screening labs. With a carefully curated genomic database established by genomic specialists, now laboratories can expand or kickstart their sequencing endeavors with confidence.

Rooted in evidence from both public and private research domains, our panel includes 390+ annotated genes across 13 different research areas (ex: immunology, metabolism, etc) enabling the detection of 87,000+ variants. We strive to offer unmatched breadth and granularity, simplifying genetic interpretation for all labs.



Stay up-to-date. Genomic discoveries arise every day - this panel is structured to reflect these findings. Our software is consistently updated with new genes and newly detected variants, whether documented by a 3rd party laboratory or by Revvity's global network of labs.* In the ever-evolving field of NGS, we help you stay current in your field of research.

Official gene panel

390+ newborn gene annotations and counting

RESEARCH AREA	GENE
Metabolism	ABCG5
Metabolism	ABCG8
Metabolism	ACAD8
Metabolism	ACAD9
Metabolism	ACADM
Metabolism	ACADS
Metabolism	ACADSB
Metabolism	ACADVL
Metabolism	ACAT1
Metabolism	AGL
Metabolism	ALDOB
Metabolism	AMT
Metabolism	ARG1
Metabolism	ARSB
Metabolism	ASL
Metabolism	ASS1
Metabolism	ATP7A
Metabolism	ATP7B
Metabolism	AUH
Metabolism	BCKDHA
Metabolism	BCKDHB
Metabolism	BTD
Metabolism	CBS
Metabolism	COQ4
Metabolism	CPS1
Metabolism	CPT1A
Metabolism	CPT2
Metabolism	CYP27A1
Metabolism	DBT
Metabolism	DHCR7
Metabolism	DLD
Metabolism	ECHS1
Metabolism	ETFA
Metabolism	ETFB
Metabolism	ETFDH
Metabolism	ETHE1
Metabolism	FAH
Metabolism	FBP1
Metabolism	G6PC1
Metabolism	GAA
Metabolism	GALC
Metabolism	GALE
Metabolism	GALK1
Metabolism	GALNS
Metabolism	GALT
Metabolism	GAMT
Metabolism	GATM
Metabolism	GBA1
Metabolism	GBE1
Metabolism	GCDH
Metabolism	GCH1
Metabolism	GLA
Metabolism	GLB1
Metabolism	GLDC
Metabolism	GLUD1
Metabolism	GSS
Metabolism	GUSB
Metabolism	GYS2
Metabolism	HADH
Metabolism	HADHA
Metabolism	HADHB
Metabolism	HCFC1
Metabolism	HGSNAT
Metabolism	HLCS
Metabolism	HMGCL
Metabolism	HMGCS2
Metabolism	HPD
Metabolism	HSD17B10
Metabolism	IDS
Metabolism	IDUA
Metabolism	IVD

RESEARCH AREA	GENE
Metabolism	LIPA
Metabolism	LMBRD1
Metabolism	LPIN1
Metabolism	MAT1A
Metabolism	MCCC1
Metabolism	MCCC2
Metabolism	MCEE
Metabolism	MLYCD
Metabolism	MMAA
Metabolism	MMAB
Metabolism	MMACHC
Metabolism	MMADHC
Metabolism	MMUT
Metabolism	MPI
Metabolism	MTHFR
Metabolism	MTR
Metabolism	MTRR
Metabolism	NADK2
Metabolism	NAGLU
Metabolism	NAGS
Metabolism	NPC1
Metabolism	OAT
Metabolism	OTC
Metabolism	OXCT1
Metabolism	PAH
Metabolism	PCBD1
Metabolism	PCCA
Metabolism	PCCB
Metabolism	PDHA1
Metabolism	PHGDH
Metabolism	PHKA2
Metabolism	PHKB
Metabolism	PHKG2
Metabolism	PIGA
Metabolism	PKLR
Metabolism	PNPO
Metabolism	POR
Metabolism	PRODH
Metabolism	PTS
Metabolism	PYGL
Metabolism	PYGM
Metabolism	QDPR
Metabolism	SGSH
Metabolism	SLC22A5
Metabolism	SLC25A13
Metabolism	SLC25A15
Metabolism	SLC25A20
Metabolism	SLC2A1
Metabolism	SLC37A4
Metabolism	SLC39A4
Metabolism	SLC7A7
Metabolism	SMPD1
Metabolism	SURF1
Metabolism	TAT
Metabolism	TCN2
Immunology	ADA
Immunology	AK2
Immunology	BLNK
Immunology	BTK
Immunology	CD247
Immunology	CD3D
Immunology	CD3E
Immunology	CD40LG
Immunology	CD79A
Immunology	CD79B
Immunology	CIITA
Immunology	CORO1A
Immunology	CYBA
Immunology	CYBB
Immunology	DCLRE1C
Immunology	DOCK8

RESEARCH AREA	GENE
Immunology	ELANE
Immunology	FOXN1
Immunology	FOXP3
Immunology	HAX1
Immunology	IGHM
Immunology	IGLL1
Immunology	IL2RG
Immunology	IL7R
Immunology	JAK3
Immunology	LIG4
Immunology	MEFV
Immunology	NCF2
Immunology	NHEJ1
Immunology	ORAI1
Immunology	PAX1
Immunology	PIK3CD
Immunology	PIK3R1
Immunology	PNP
Immunology	PRF1
Immunology	PRKDC
Immunology	PTPRC
Immunology	RAC2
Immunology	RAG1
Immunology	RAG2
Immunology	RFX5
Immunology	RFXANK
Immunology	RFXAP
Immunology	SLC39A7
Immunology	STAT3
Immunology	STIM1
Immunology	TCF3
Immunology	TOP2B
Immunology	UNC13D
Immunology	WAS
Immunology	XIAP
Immunology	ZAP70
Cardiovascular	ACTA2
Cardiovascular	ACTC1
Cardiovascular	AOB
Cardiovascular	BAG3
Cardiovascular	CALM1
Cardiovascular	CALM2
Cardiovascular	CALM3
Cardiovascular	CASQ2
Cardiovascular	CHD7
Cardiovascular	COL3A1
Cardiovascular	DES
Cardiovascular	DSC2
Cardiovascular	DSG2
Cardiovascular	DSP
Cardiovascular	FBN1
Cardiovascular	KCNE1
Cardiovascular	KCNH2
Cardiovascular	KCNQ1
Cardiovascular	LDLR
Cardiovascular	LMNA
Cardiovascular	LPL
Cardiovascular	MYBPC3
Cardiovascular	MYH11
Cardiovascular	MYH7
Cardiovascular	MYL2
Cardiovascular	MYL3
Cardiovascular	PCSK9
Cardiovascular	PKP2
Cardiovascular	PRKAG2
Cardiovascular	PTPN11
Cardiovascular	RBM20
Cardiovascular	RYR2
Cardiovascular	SCN5A
Cardiovascular	SMAD3
Cardiovascular	SMAD4

RESEARCH AREA	GENE
Cardiovascular	SOS1
Cardiovascular	TAFAZZIN
Cardiovascular	TBX1
Cardiovascular	TBX2
Cardiovascular	TGFBR1
Cardiovascular	TGFBR2
Cardiovascular	TMEM43
Cardiovascular	TNNC1
Cardiovascular	TNNI3
Cardiovascular	TNNT2
Cardiovascular	TPM1
Cardiovascular	TRDN
Cardiovascular	TTN
Cardiovascular	ZIC3
Endocrinology	ABCC8
Endocrinology	ALPL
Endocrinology	ANOS1
Endocrinology	AQP2
Endocrinology	AVPR2
Endocrinology	CASR
Endocrinology	CYP11B1
Endocrinology	CYP11B2
Endocrinology	CYP17A1
Endocrinology	CYP27B1
Endocrinology	DUOX2
Endocrinology	DUOXA2
Endocrinology	GH1
Endocrinology	GHR
Endocrinology	GHRHR
Endocrinology	HNF1A
Endocrinology	HNF4A
Endocrinology	HSD3B2
Endocrinology	IGSF1
Endocrinology	INS
Endocrinology	INSR
Endocrinology	IYD
Endocrinology	KCNJ11
Endocrinology	LHX3
Endocrinology	MC2R
Endocrinology	NROB1
Endocrinology	PAX8
Endocrinology	PHEX
Endocrinology	POU1F1
Endocrinology	PROKR2
Endocrinology	PROP1
Endocrinology	SCNN1A
Endocrinology	SCNN1B
Endocrinology	SLC34A3
Endocrinology	SLC5A5
Endocrinology	SRY
Endocrinology	STAR
Endocrinology	TCIRG1
Endocrinology	TG
Endocrinology	THRA
Endocrinology	TPO
Endocrinology	TRHR
Endocrinology	TSHB
Endocrinology	TSHR
Endocrinology	VDR
Neurology	ABCD1
Neurology	ALDH7A1
Neurology	ARSA
Neurology	ATM
Neurology	BCAP31
Neurology	CACNA1S
Neurology	CLCN1
Neurology	CLN3
Neurology	CRLF1
Neurology	DMD
Neurology	FLAD1
Neurology	FOLR1

RESEARCH AREA	GENE
Neurology	KCNQ2
Neurology	NF1
Neurology	NTRK1
Neurology	PRRT2
Neurology	RYR1
Neurology	SCN1A
Neurology	SCN2A
Neurology	SCN8A
Neurology	SLC52A2
Neurology	SLC52A3
Neurology	SPR
Neurology	TH
Neurology	TSC1
Neurology	TSC2
Neurology	TTPA
Hematology	ACVRL1
Hematology	ADAMTS13
Hematology	ANK1
Hematology	DKC1
Hematology	ENG
Hematology	EPB42
Hematology	F9
Hematology	FANCA
Hematology	G6PD
Hematology	HBB
Hematology	HBD
Hematology	HPS1
Hematology	HPS4
Hematology	LYST
Hematology	MPL
Hematology	RPL11
Hematology	RPS19
Hematology	RPS26
Hematology	SBDS
Hematology	SH2D1A
Hematology	SPTB
Hematology	TERT
Hematology	VWF
Oncology	APC
Oncology	BMPR1A
Oncology	ERCC2
Oncology	ERCC5
Oncology	MAX
Oncology	MEN1
Oncology	NF2
Oncology	PTCH1
Oncology	PTEN
Oncology	RB1
Oncology	RET
Oncology	SDHAF2
Oncology	SDHB
Oncology	SDHC
Oncology	SDHD
Oncology	STK11
Oncology	TMEM127
Oncology	TP53
Oncology	VHL
Oncology	WT1
Oncology	XPA
Oncology	XPC
ENT/Dental	CDH23
ENT/Dental	GJB2
ENT/Dental	GJB6
ENT/Dental	MYO15A
ENT/Dental	PAX3
ENT/Dental	SLC26A4
ENT/Dental	SOX10
ENT/Dental	TECTA
ENT/Dental	TMIE
ENT/Dental	TMPRSS3
ENT/Dental	TPRN

RESEARCH AREA	GENE
ENT/Dental	TRIOBP
ENT/Dental	USH1C
ENT/Dental	USH1G
ENT/Dental	USH2A
ENT/Dental	WHRN
Nephrology	AGXT
Nephrology	CFH
Nephrology	COL4A3
Nephrology	COL4A4
Nephrology	COL4A5
Nephrology	CTNS
Nephrology	GRHRP
Nephrology	HOGA1
Nephrology	PKD2
Nephrology	PKHD1
Nephrology	SLC12A3
Nephrology	SLC4A1
Gastroenterology	ABCB11
Gastroenterology	ABCB4
Gastroenterology	AKR1D1
Gastroenterology	ATP8B1
Gastroenterology	HSD3B7
Gastroenterology	IL10RA
Gastroenterology	JAG1
Gastroenterology	MTPP
Gastroenterology	SLCO1B1
Gastroenterology	TRMU
Gastroenterology	UGT1A1
Ophthalmology	CYP1B1
Ophthalmology	GPR143
Ophthalmology	OPA1
Ophthalmology	RPE65
Ophthalmology	SLC45A2
Orthopedics	ANKH
Orthopedics	COL1A1
Orthopedics	COL1A2
Orthopedics	RMRP
Dermatology	KRT14
Dermatology	KRT5
Pulmonology	CFTR

For research use only. Not for use in diagnostic procedures.



For research use only. Not for
use in diagnostic procedures.

Kickstart your genomic discoveries today.

Now you can access Revvity's expansive gene panel through our comprehensive NGS software. Beyond interpretation of DNA data, this solution also includes LIMS and DNA analysis capabilities, backed with the latest data security protocols. See how you can amplify your research alongside the global leaders in newborn screening today.

[Learn More](#)

revvity