



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 100,000+ variants. We strive to offer deeper insights while simplifying genomic analysis for all labs.

Stay up-to-date

Genomic discoveries arise every day – this panel is structured to reflect these findings. Consistently updated with new genes and newly detected variants, whether documented by a 3rd party laboratory or by Revvity’s global network of labs,* we help you stay current in your field of research.

*Revvity Omics lab locations: USA, UK, India, Sweden, China

Official gene panel: 390+ newborn gene annotations and counting

Metabolism

ABCG5	ATP7A	ETFA	GBE1	HMGCL	MMAB	PCBD1	QDPR
ABCG8	ATP7B	ETFB	GCDH	HMGCS2	MMACHC	PCCA	SGSH
ACAD8	AUH	ETFDH	GCH1	HPD	MMADHC	PCCB	SLC22A5
ACAD9	BCKDHA	ETHE1	GLA	HSD17B10	MMUT	PDHA1	SLC25A13
ACADM	BCKDHB	FAH	GLB1	IDS	MPI	PHGDH	SLC25A15
ACADS	BTD	FBP1	GLDC	IDUA	MTHFR	PHKA2	SLC25A20
ACADSB	CBS	G6PC1	GLUD1	IVD	MTR	PHKB	SLC2A1
ACADVL	COQ4	GAA	GSS	LIPA	MTRR	PHKG2	SLC37A4
ACAT1	CPS1	GALC	GUSB	LMBRD1	NADK2	PIGA	SLC39A4
AGL	CPT1A	GALE	GYS2	LPIN1	NAGLU	PKLR	SLC7A7
ALDOB	CPT2	GALK1	HADH	MAT1A	NAGS	PNPO	SMPD1
AMT	CYP27A1	GALNS	HADHA	MCCC1	NPC1	POR	SURF1
ARG1	DBT	GALT	HADHB	MCCC2	OAT	PRODH	TAT
ARSB	DHCR7	GAMT	HCFC1	MCEE	OTC	PTS	TCN2
ASL	DLD	GATM	HGSNAT	MLYCD	OXCT1	PYGL	
ASS1	ECHS1	GBA1*	HLCS	MMAA	PAH	PYGM	

Immunology

ADA	CD40LG	DCLRE1C	IGLL1	NHEJ1	PRKDC	RFXAP	WAS
AK2	CD79A	DOCK8	IL2RG	ORAI1	PTPRC	SLC39A7	XIAP
BLNK	CD79B	ELANE	IL7R	PAX1	RAC2	STAT3	ZAP70
BTK	CIITA	FOXN1	JAK3	PIK3CD	RAG1	STIM1	
CD247	CORO1A	FOXP3	LIG4	PIK3R1	RAG2	TCF3	
CD3D	CYBA	HAX1	MEFV	PNP	RFX5	TOP2B	
CD3E	CYBB	IGHM	NCF2	PRF1	RFXANK	UNC13D	

Cardiovascular

ACTA2	CASQ2	FBN1	MYBPC3	PRKAG2	SOS1	TNNC1
ACTC1	CHD7	KCNE1	MYH11	PTPN11	TAFAZZIN	TNNI3
APOB	COL3A1	KCNH2	MYH7	RBM20	TBX1	TNNT2
BAG3	DES	KCNQ1	MYL2	RYR2	TBX2	TPM1
CALM1	DSC2	LDLR	MYL3	SCN5A	TGFBR1	TRDN
CALM2	DSG2	LMNA	PCSK9	SMAD3	TGFBR2	TTN
CALM3	DSP	LPL	PKP2	SMAD4	TMEM43	ZIC3

Endocrinology

ABCC8	CYP11B1	GH1	IGSF1	MC2R	PROP1	STAR	TSHB
ALPL	CYP11B2	GHR	INS	NR0B1	SCNN1A	TCIRG1	TSHR
ANOS1	CYP17A1	GHRHR	INSR	PAX8	SCNN1B	TG	VDR
AQP2	CYP27B1	HNF1A	IYD	PHEX	SLC34A3	THRA	
AVPR2	DUOX2	HNF4A	KCNJ11	POU1F1	SLC5A5	TPO	
CASR	DUOXA2	HSD3B2	LHX3	PROKR2	SRY	TRHR	

Neurology

ABCD1*	BCAP31	CRLF1	KCNQ2	RYR1	SLC52A2	TSC1
ALDH7A1	CACNA1S	DMD	NF1	SCN1A	SLC52A3	TSC2
ARSA	CLCN1	FLAD1	NTRK1	SCN2A	SPR	TTPA
ATM	CLN3	FOLR1	PRRT2	SCN8A	TH	

Hematology

ACVRL1	DKC1	F9	HBB	HPS4	RPL11	SBDS	TERT
ADAMTS13	ENG	FANCA	HBD	LYST	RPS19	SH2D1A	VWF
ANK1	EPB42	G6PD	HPS1	MPL	RPS26	SPTB	

Oncology

APC	ERCC5	NF2	RB1	SDHB	STK11	VHL	XPC
BMPR1A	MAX	PTCH1	RET	SDHC	TMEM127	WT1	
ERCC2	MEN1	PTEN	SDHAF2	SDHD	TP53	XPA	

ENT/Dental

CDH23	GJB6	PAX3	SOX10	TMIE	TPRN	USH1C	USH2A
GJB2	MYO15A	SLC26A4	TECTA	TMPRSS3	TRIOBP	USH1G	WHRN

Nephrology

AGXT	COL4A4	GRHPR	PKHD1
CFH	COL4A5	HOGA1	SLC12A3
COL4A3	CTNS	PKD2	SLC4A1

Gastroenterology

ABCB11	HSD3B7	SLCO1B1
ABCB4	IL10RA	TRMU
AKR1D1	JAG1	UGT1A1
ATP8B1	MTTP	

Orthopedics

ANKH
COL1A1
COL1A2
RMRP

Ophthalmology

CYP1B1	OPA1	SLC45A2
GPR143	RPE65	

Dermatology

KRT14
KRT5

Pulmonology

CFTR

* Long-range PCR-kit (sold separately) is required for the library preparation step of highlighted genes.

* Long-range PCR-kit (sold separately) is required for the library preparation step of highlighted genes.

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.



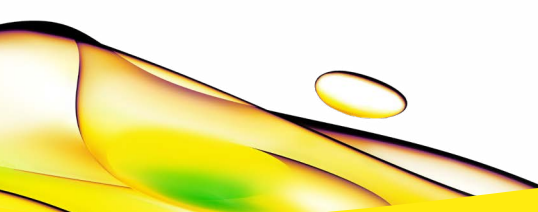
[Learn more](#)

Scan or click
for more information



[Contact us](#)

Scan or click
to contact us.



revvity

www.revvity.com

Copyright ©2025, Revvity, Inc. All rights reserved.
Revvity is a trademark of Revvity, Inc. All other trademarks are the property of their respective owners.

2133350

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