



# superRCA On-demand Assay Catalog

These assays are available and will be prepared after order. The assay name should be added as a comment in quotes and orders.

Assay name	Amino Acid Change	cDNA Mutation	DNA Input
AKT1_Q545L	p.Q545L		66ng
ARID1A_S2264*	p.S2264*		66ng
ASXL1 F808LfsX10	F808LfsX10		660ng
ASXL1 p.R693Ter	p.R693Ter	c.2077C>T	660ng
ASXL1_C687fs30	p.C687fs30		660ng
ASXL1_L764fs8	p.L764fs8		660ng
ASXL1_S1146x	p.S1146x		660ng
ATM_p.Arg2832Cys	p.Arg2832Cys		66ng
ATP1A1 F100_L104del	p.F100_L104del	c.299_313del	66ng
ATP1A1 L104R	p.L104R	c.311T>G	66ng
ATP2B3 L425_V426del	p.L425_V426del	c.1272_1277del	66ng
ATP2B3 L425_V426del	p.L425_V426del	c.1273_1278del	66ng
BCOR p.Arg1183Ter	p.Arg1183Ter		660ng
BCOR p.Tyr1548Ter	p.Tyr1548Ter		660ng
BRAF V600E	p.V600E		
BRAF V600E	p.V600E		
BRAF V600K	p.V600K		
BRAF V600K	p.V600K		
BRAF_G509A	p.G509A		66ng
BRAF_p.Asn621Ile	p.Asn621Ile		66ng
BRAF_p.Leu637Gln	p.Leu637Gln		66ng
BRCA2_Q969*	p.Q969*		66ng
CACNA1D p.F747V/c.2239T>C	p.F747V		66ng
CACNA1D p.F747V/c.2239T>G	p.F747V		66ng
CACNA1D p.G403R	p.G403R		66ng
CARD11_p.Met360Val	p.Met360Val		66ng
CBL C381Y	p.C381Y		660ng
CBL L380P	p.L380P		660ng
CBL p.Trp408Cys	p.Trp408Cys		660ng
CDH20_S99	N/A		66ng
CDK12_G871V	p.G871V		66ng
CTCF_D443N	p.D443N		66ng
DDX41 p.Arg525His	p.Arg525His		660ng
DNMT3A_2597+1G/del			660ng
DNMT3A_Q816R	p.Q816R		660ng
DNMT3A_R379C	p.R379C		660ng
EFGR T790M	p.T790M		
EGFR 2235-2249del			
EGFR 2235-2250del			
EGFR 2236-2250del			
EGFR 2239-2248delinsC			



Assay name	Amino Acid Change	cDNA Mutation	DNA Input
EGFR 2240-2254del			
EGFR C797SA	p.797SA		
EGFR C797SC	p.797SC		
EGFR G719S	p.G719S		
EGFR L858R	p.L858R		
EGFR L861Q	p.L861Q		
EGFR_P753L	p.P753L		
EP300_Q102R	p.Q102R		66ng
ERCC6_E1351Q	p.E1351Q		66ng
ESR1_D538G	D538G		66ng
ESR1_E380Q	E380Q		66ng
ESR1_L536H	L536H		66ng
ESR1_L536P	L536P		66ng
ESR1_L536R	L536R		66ng
ESR1_S463P	S463P		66ng
ESR1_V422del	V422del		66ng
ESR1_Y537C	Y537C		66ng
ESR1_Y537N	Y537N		66ng
ESR1_Y537S	Y537S		66ng
ETV6 L348F	p.L348F		660ng
ETV6 p.Leu398Pro	p.Leu398Pro		660ng
EZH2_R684C	p.R684C		660ng
EZH2_Y733Lfs5			660ng
FAT1_p.Asp1956Asn	p.Asp1956Asn		66ng
FLT3 p.Asp835His	p.Asp835His		660ng
FMC63CAR/GAPDH			660ng
GATA2 G366D367dup	G366D367dup		660ng
HRAS_M111fs	p.M111fs		66ng
IDH1_R132C	p.R132C		660ng
IDH1_R132G	p.R132G		660ng
IDH1_R132H	p.R132H		660ng
IDH1_R132L	p.R132L		660ng
IDH1_R132S	p.R132S		660ng
IDH2_R140L	p.R140L		660ng
IDH2_R140Q	p.R140Q		660ng
IDH2_R140W	p.R140W		660ng
IDH2_R172K	p.R172K		660ng
IDH2_R172M	p.R172M		660ng
IGLL5_p.Gly6Asp	p.Gly6Asp		66ng
JAK2_V617F	p.V617F		660ng
JAK3 p.R657N	p.R657N		660ng
KCNJ5 G151R	p.G151R		66ng
KCNJ5 L168R	p.L168R		66ng
KIT_D816V	p.D816V		660ng
KIT_S285C	p.S285C		66ng
KRAS A146V	p.A146V		
KRAS A59T	p.A59T		
KRAS A59T	p.A59T		
KRAS G12A	p.G12A		
KRAS G12A	p.G12A		
KRAS G12C	p.G12C		
KRAS G12C	p.G12C		
KRAS G12D	p.G12D		
KRAS G12D	p.G12D		
KRAS G12E	p.G12E		
KRAS G12E	p.G12E		
KRAS G12R	p.G12R		
KRAS G12R	p.G12R		
KRAS G12S	p.G12S		
KRAS G12S	p.G12S		
KRAS G12V	p.G12V		

Assay name	Amino Acid Change	cDNA Mutation	DNA Input
KRAS G12V	p.G12V		
KRAS G13A	p.G13A		
KRAS G13C	p.G13C		
KRAS G13C	p.G13C		
KRAS G13D	p.G13D		
KRAS G13D	p.G13D		
KRAS G13R	p.G13R		
KRAS G13R	p.G13R		
KRAS G13S	p.G13S		
KRAS G13V	p.G13V		
KRAS K147Q	p.K147Q		
KRAS Q61Hc	p.Q61Hc		
KRAS Q61Hc	p.Q61Hc		
KRAS Q61Ht	p.Q61Ht		
KRAS Q61Ht	p.Q61Ht		
KRAS Q61L	p.Q61L		
KRAS Q61L	p.Q61L		
KRAS_G12C	p.G12C		66ng
KRASA146T	p.A146T		
MAP2K4_R315*	p.R315*		66ng
MBD1_E757E	p.E757E		66ng
MET_D1010Y	p.D1010Y		66ng
MET_S889F	p.S889F		66ng
MIR6871_n.*4768A>C	n.*4768A>C		66ng
MYD88_L265P	L265P		660ng
NF1_Y489C	p.Y489C		66ng
NFE2L2_Q26_D27insVLWRQ	N/A		66ng
NPM1_MutA (TCTG/dup)	p.W288Cfs*12.		660ng
NPM1_MutB (CATG)	p.W288Cfs*12		660ng
NPM1_MutD (CCTG)	p.W288Cfs*12		660ng
NRAS A59T	p.A59T		66ng
NRAS G12A	p.G12A		66ng
NRAS G12C	p.G12C		66ng
NRAS G12D	p.G12D		66ng
NRAS G12R	p.G12R		66ng
NRAS G12S	p.G12S		66ng
NRAS G12V	p.G12V		66ng
NRAS G13D	p.G13D		66ng
NRAS G13R	p.G13R		66ng
NRAS G13V	p.G13V		66ng
NRAS Q61Hc	p.Q61Hc		66ng
NRAS Q61Ht	p.Q61Ht		66ng
NRAS Q61K	p.Q61K		66ng
NRAS Q61L	p.Q61L		66ng
NRAS Q61P	p.Q61P		66ng
NRAS Q61R	p.Q61R		66ng
NSD2_p.Thr1150Ala	p.Thr1150Ala		66ng
PARP4_c.3367-4A>C	N/A		66ng
PIK3C2G_I169M	p.I169M		66ng
PIK3CA E542K	p.E542K		66ng
PIK3CA E545G	p.E545G		66ng
PIK3CA E545K	p.E545K		66ng
PIK3CA H1047L	p.H1047L		66ng
PIK3CA H1047R	p.H1047R		66ng
PIK3CA H1047Y	p.H1047Y		66ng
PIK3CA I112F	p.I112F		66ng
PIK3CA K111E	p.K111E		66ng
PIK3CA M1043I	p.M1043I		66ng
PIK3CA Q546K	p.Q546K		66ng
PIK3CA Q546R	p.Q546R		66ng
PTEN_Q214*	p.Q214*		66ng

Assay name	Amino Acid Change	cDNA Mutation	DNA Input
PTEN_S229*	p.S229*		66ng
PTPN11 p.E58Y		c.172A>T	660ng
PTPN11 p.E59Y	p.E59Y	c.205G>A	660ng
PTPN11_S502P	p.S502P		660ng
RB1_I103fs	p.I103fs		66ng
RET_G1071E	p.G1071E		66ng
ROS1_Q379R	p.Q379R		66ng
RUNX1 p.Arg107Cys	p.Arg107Cys		660ng
RUNX1 p.Arg166Gly	p.Arg166Gly		660ng
RUNX1_Ser322X	p.Ser322X		660ng
RUNX1_Val164SerfsX48	p.Val164SerfsX48		660ng
SETBP1 p.D868N	p.D868N	c.2602G>A	660ng
SETBP1_Asp868Asn	p.Asp868Asn		660ng
SF3B1_L700E	p.L700E		660ng
STAG2 R590*	p.R590*		660ng
STK11_Q100E	p.Q100E		66ng
TAF1_R1238Q	p.R1238Q		66ng
TERT_C228T	p.C228T		
TERT_C250T	p.C250T		
TET2_C1709fs10	p.C1709fs10		660ng
TET2_F1104fs*3	p.F1104fs*3		660ng
TET2_R1261C	p.R1261C		660ng
TET2_T1895fs*11	p.T1895fs*11		660ng
TP53_260del*	N/A		66ng
TP53_c.673-2A>T	N/A		66ng
TP53_c.783-2A>C	N/A		66ng
TP53_C238Y	p.C238Y		66ng
TP53_E294*	p.E294*		66ng
TP53_E339*	p.E339*		66ng
TP53_H179R	p.H179R		66ng
TP53_p.Phe113Cys	p.Phe113Cys		66ng
TP53_V173L	p.V173L		66ng
TP53_V272E	p.V272E		66ng
U2AF1_Q157P	p.Q157P		

For Research Use Only. Not for use in diagnostic procedures.  
superRCA is a trademark or registered trademark of Rarity Bioscience AB.

© 2025 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries. All other trademarks are the property of their respective owners.

For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at [beckman.com](https://www.beckman.com)  
2025-GBL-EN-108190-v1

