



TREVIGEN®

Tools to Study Genomic Instability and Genotoxic Stress DNA RepairGene Knockdown Cell Lines Base Excision Repair (BER)

Trevigen's Base Excision Repair Knockdown Cell Lines are target specific LN428 (glioblastoma) shRNA lentivirus transduced cells. They are rigorously qualified and mycoplasma free. The percent knockdown levels are evaluated by RT-PCR. Lentiviruses are maintained by puromycin selection.

Ordering information*

Target	Product Name (Knockdown cell line)	Catalog Number	%KD by RT-PCR	# of cells per vial	Price
APE1	KD-BER-LN428-APE1	5517-001-01 ◊	90%	1x10 ⁶	\$895
APE2	KD-BER-LN428-APE2	5518-001-01	80%	1x10 ⁶	\$895
BRCA1	KD-HR-LN428-BRCA1**	5502-001-01	83%	1x10 ⁶	\$895
XRCC1	KD-BER-LN428-XRCC1	5516-001-01	81%	1x10 ⁶	\$895
MBD4	KD-BER-LN428-MBD4	5506-001-01	72%	1x10 ⁶	\$895
MPG	KD-BER-LN428-MPG	5511-001-01	98%	1x10 ⁶	\$895
MutYH	KD-BER-LN428-MutYH	5512-001-01	87%	1x10 ⁶	\$895
NEIL1	KD-BER-LN428-NEIL1	5513-001-01	92%	1x10 ⁶	\$895
NEIL2	KD-BER-LN428-NEIL2	5507-001-01	86%	1x10 ⁶	\$895
NEIL3	KD-BER-LN428-NEIL3	5508-001-01	95%	1x10 ⁶	\$895
NTHL1	KD-BER-LN428-NTHL1	5505-001-01 ◊	91%	1x10 ⁶	\$895
OGG1	KD-BER-LN428-OGG1	5504-001-01	63%	1x10 ⁶	\$895
PARG	KD-BER-LN428-PARG	5501-001-01	84%	1x10 ⁶	\$895
PARP1	KD-BER-LN428-PARP1	5500-001-01	72%	1x10 ⁶	\$895
PARP2	KD-BER-LN428-PARP2	5514-001-01	83%	1x10 ⁶	\$895
PARP3	KD-BER-LN428-PARP3	5515-001-01	70%	1x10 ⁶	\$895
SMUG1	KD-BER-LN428-SMUG1	5510-001-01	63%	1x10 ⁶	\$895
TDG	KD-BER-LN428-TDG	5519-001-01	74%	1x10 ⁶	\$895
UNG	KD-BER-LN428-UNG	5509-001-01	87%	1x10 ⁶	\$895
Control	KD-BER-LN428-Control	5503-001-01	N/A	1x10 ⁶	\$475

* Cell lines are provided under a Limited License MTA. Please inquire at info@trevigen.com.

** BRCA1 participates in the homologous recombination pathway and shows synthetic lethality with PARP1.

◊ Cell lines available soon.

Successful and consistent results are dependent upon the quality and degree of characterization of the cells under investigation. Highly passaged cells may undergo both genotypic and phenotypic changes that render them an inadequate *in vitro* model for specific investigations. We recommend for all studies that highly qualified low passage number cells are used to ensure reliable and reproducible results.

Validation of BER Knockdown cell lines

ANGIOGENESIS • DNA DAMAGE & REPAIR • APOPTOSIS

Trevigen, Inc.
8405 Helgerman Court
Gaithersburg, Maryland 20877
1-800-873-8443



TREVIGEN®

Tools to Study Genomic Instability and Genotoxic Stress

DNA RepairGene Knockdown Cell Lines Base Excision Repair (BER)

All Trevigen knockdown cell lines are evaluated by RT-PCR and Western Blots when reliable immunoreagents are available. Additional functional assays are performed when feasible.

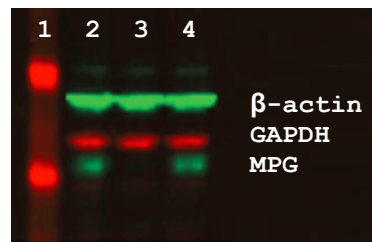
ANGIOGENESIS • DNA DAMAGE & REPAIR • APOPTOSIS

Enzymatic Assay for MPG activity



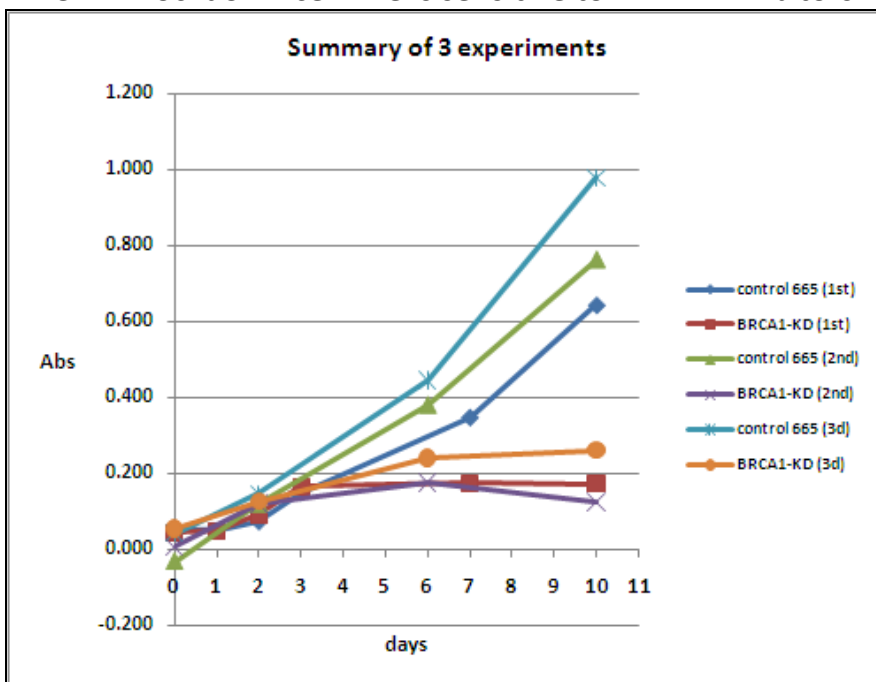
Lane 1: Buffer alone
Lane 2: Control extract
Lane 3: MPG knockdown extract

Western Blot of MPG protein



Lane 1: Molecular weight markers
Lane 2: Scramble RNA Control
Lane 3: MPG knockdown extract
Lane 4: Control cell line

BRCA1 Knockdown cell line is sensitive to PARP1 inhibitors



©2011 Trevigen is a registered trademark of Trevigen, Inc.

Trevigen, Inc.
8405 Helgerman Court
Gaithersburg, Maryland 20877
1-800-873-8443