



## LIST OF KNOCKDOWN CELL LINES

### Base Excision Repair

Catalog #	Product Name	# of Cells Per Vial	Size	Price	%KD by RT-PCR
5500-001-01	PARP1 Knock Down Cell Line (MTA Required prior to shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	72%
5504-001-01	OGG1 Knock Down Cell Line (MTA Required prior to shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	63%
5505-001-01	NTHL1 Knock Down Cell Line (MTA Required prior to shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	91%
5507-001-01	NEIL2 Knock Down Cell Line (MTA Required prior to shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	86%
5508-001-01	NEIL3 Knock Down Cell Line (MTA Required prior to shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	95%
5509-001-01	UNG Knock Down Cell Line (MTA Required prior to shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	87%
5510-001-01	SMUG1 Knock Down Cell Line (MTA Required prior to shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	63%
5511-001-01	MPG Knock Down Cell Line (MTA Required prior to shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	98%
5512-001-01	MutYH Knock Down Cell Line (MTA Required prior to shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	87%
5513-001-01	Neil1 Knock Down Cell Line (MTA Required prior to shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	92%
5514-001-01	PARP2 Knock Down Cell Line (MTA Required prior to shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	83%

	to shipment)				
5515-001-01	PARP3 Knock Down Cell Line (MTA Required prior to shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	70%
5516-001-01	XRCC1 Knock Down Cell Line (MTA Required prior to shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	81%
5517-001-01	APE1 Knock Down Cell Line (MTA Required prior to shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	90%
5518-001-01	APE2 Knock Down Cell Line (MTA Required prior to shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	80%
5519-001-01	TDG Knock Down Cell Line (MTA Required prior to shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	74%
5506-001-01	MBD4 Knock Down Cell Line (MTA required before shipment)	1x10	1 vial	\$895.00	72%
54999-001-01	LIG3 Knock Down Cell Line (MTA required prior to shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	88%

## Mismatch Repair

Catalog #	Product Name	# of Cells Per Vial	Size	Price	%KD by RT-PCR
55020-001-01	MSH2 Knock down cell line (MTA required)	1x10 <sup>6</sup>	1 vial	\$895.00	70%
55027-001-01	PMS2 Knock Down Cell Line (MTA Required prior to shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	84%
55028-001-01	MLH3 Knock Down Cell Line (MTA required before shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	71%
55021-001-01	MSH3 Knock Down Cell Line (MTA required before shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	83%
55026-001-01	MLH1 Knock Down Cell Lines (MTA required to shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	80%
55024-001-01	MSH5 Knock Down Cell Line (MTA required to shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	92%

55022-001-01	MSH6 Knock Down Cell Line (MTA required to shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	93%
--------------	--	-------------------	--------	----------	-----

## Non-Homologous End Joining

Catalog #	Product Name	# of Cells Per Vial	Size	Price	%KD by RT-PCR
55041-001-01	SIRT5 Knock down cell line (MTA required)	1x10 <sup>6</sup>	1 vial	\$895.00	83%
55037-001-01	SIRT1 Knock Down Cell Line (MTA required prior to shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	95%
55036-001-01	PRKDC Knock Down Cell Line (MTA required prior to shipping)	1x10 <sup>6</sup>	1 vial	\$895.00	94%
55038-001-01	SIRT2 Knock Down Cell Line (MTA required prior to shipping)	1x10 <sup>6</sup>	1 vial	\$895.00	98%
55034-001-01	XRCC6 (Ku70) Knock Down Cell Line (MTA required before shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	92%
55035-001-01	XRCC5 (Ku80) Knock Down Cell Line (MTA required before shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	85%
55039-001-01	SIRT3 Knock Down Cell Line (MTA required)	1 x 10 <sup>6</sup>	1 vial	\$895.00	86%

## Knockdown Control

Catalog #	Product Name	# of Cells Per Vial	Size	Price	%KD by RT-PCR
5503-001-01	Knock Down Control Cell Line (MTA Required prior to shipment)	1x10 <sup>6</sup>	1 vial	\$475.00	N/A

## Homologous Recombination

Catalog #	Product Name	# of Cells Per Vial	Size	Price	%KD by RT-PCR
-----------	--------------	---------------------	------	-------	---------------

55056-001-01	RAD54B Knock down cell line (MTA required prior to shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	80%
55062-001-01	NBS1 Knock Down Cell Line (MTA required prior to shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	73%
55066-001-01	RAD21 Knock Down Cell Line (MTA required prior to shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	93%
55053-001-01	XRCC3 Knock Down Cell Line (MTA required prior to shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	93%
55065-001-01	SMC6L1 Knock Down Cell Line (MTA required prior to shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	91%
55059-001-01	SHFM1 Knock Down Cell Line (MTA required before shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	84%
5502-001-01	BRCA1 Knock Down Cell Line (MTA Required prior to shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	83%
55049-001-01	RAD51C Knock down cell line (MTA required before shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	92%
55060-001-01	RAD50 Knock Down Cell Line (MTA required prior to shipment)	1x10 <sup>6</sup>	1 vial	\$895.00	81%

### Trevigen, Inc.

8405 Helgerman Court, Gaithersburg, MD 20877 USA

Phone: 1-800-TREVIGEN (1-800-873-8443) • 301-216-2800 • Fax: 301-560-4973 • Email: [info@trevigen.com](mailto:info@trevigen.com) • <http://www.trevigen.com>