

Product Information

MISSION® shRNA Human Gene Family Sets , DNA Format

Catalog Numbers **SH0121, SH0221, SH0421, SH0521, SH0721, SH0821, SH1021, SH1121, SH1321, SH1821, SH1921, SH2121, SH2221, SH2321, SH2421, SH2521, SH2621, SH2721, SH2821, SH2921, SH3021**

Storage Temperature –20 °C

TECHNICAL BULLETIN

Product Description

Small interfering RNAs (siRNAs) generated from short hairpin RNAs (shRNAs) are a powerful way to mediate gene specific RNA interference (RNAi) for extended periods of time in mammalian cells. The MISSION product line is a viral-vector-based RNAi library against annotated mouse and human genes. MISSION shRNAs are expressed intracellularly after transduction with amphotropic lentivirus particles, allowing screening in a wide range of mammalian cell lines. In these cell lines, MISSION shRNA clones permit rapid, cost efficient loss-of-function and genetic interaction screens. We have included a table of reviews for each gene family set.

The MISSION shRNA Human Gene Family Sets, DNA Format, contains shRNA clones that allows for high throughput loss-of-function and genetic interaction screens. Each MISSION shRNA clone is constructed within the lentivirus plasmid vector pLKO.1-Puro.¹ The pLKO.1-Puro vector contains the ampicillin and puromycin antibiotic resistance genes for selection of inserts in bacterial or mammalian cells respectively. The set consists of sequence-verified shRNA lentiviral plasmid DNA. Each target set consists of 3 or more constructs that have been designed against each target gene using a proprietary algorithm. Therefore, a range of gene silencing efficiencies, with at least one construct from each gene set being >70%, can be expected when using these clones. This allows one to examine the effect of loss of gene function over a large series of gene knockdown efficiencies. Each shRNA construct has been cloned and sequence verified to ensure a match to the target gene.

RNAi knockdown be achieved either with the plasmid DNA or the MISSION lentiviral delivery system. Target cell lines may be transfected with the purified plasmid for transient or stable gene silencing (puromycin selection). In addition, self-inactivating replication

incompetent viral particles can be produced in packaging cells (HEK293T) by co-transfection with compatible packaging plasmids.²⁻³ Unlike murine-based MMLV or MSCV retroviral systems, lentiviral-based particles permit efficient infection and genomic integration of the specific shRNA construct into differentiated and non-dividing cells, such as neurons and dendritic cells, overcoming low transfection and integration difficulties when using these cell lines.

Please see the **Cell Type Table** for those cell types that have been successfully infected by pLKO.1-puro based shRNA constructs.

Components/Reagents

The individual constructs are provided as 40 µL frozen stocks containing an average of 2 µg of plasmid DNA in Tris-EDTA (TE) buffer with amounts of DNA ranging from 400 ng to 4 µg per well. Sets are provided in 96-well plates with a one dimensional barcode label on each plate and a CD containing plate map positions.

The hairpin sequence and other unique clone information may be obtained by searching the MISSION search database at: www.sigma.com/yfq using RefSeq accession numbers, e.g. NM_027088, unique clone identification numbers, e.g. NM_027088.1-989s1c1, or TRC numbers, e.g. TRCN0000030720.

Precautions and Disclaimer

These products are for R&D use only, not for drug, household, or other uses. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

Storage/Stability

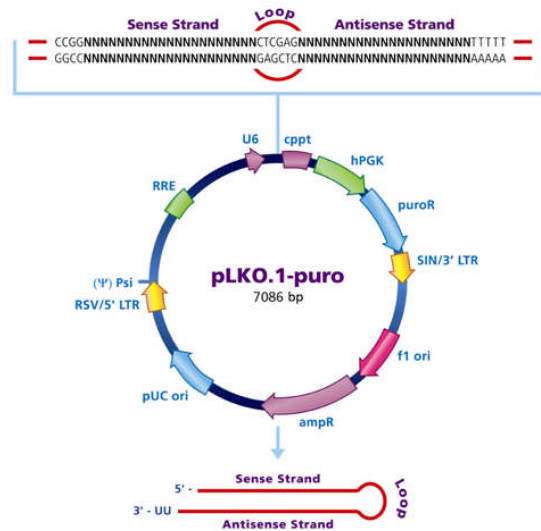
All components are stable for at least one year after receipt when stored at –20 °C.

Catalog Number	Gene Family Set	Gene Count *	Clone Count *	Average Number Clones/Gene *
SH1921	Apoptosis Pathway	440	2882	6.6
SH2921	B-Cell Activation	99	563	5.7
SH2221	Cell Adhesion Genes	366	2013	5.5
SH0821	Cytokine and Chemokine	106	525	5.0
SH1321	Cytokine and Chemokine Receptors	94	540	5.7
SH2321	Cytoskeleton Genes	273	1534	5.6
SH3021	Epigenetic Regulators	10	55	5.5
SH1821	DNA Repair Pathway	116	658	5.7
SH0721	Ubiquitin Hydrolases (DUBS)	124	612	4.9
SH2521	Extracellular Matrix Genes	330	1750	5.3
SH0221	G-Protein Coupled Receptors (GPCRs)	444	2761	6.2
SH2621	Helicase	133	676	5.1
SH1021	Ion Channel	276	1357	4.9
SH2721	JAK-STAT Pathway	189	1128	6.0
SH0121	Kinases, complete	513	5141	10.0
SH1121	Nuclear Hormone Receptors	46	218	4.7
SH2421	p53 Pathway	239	1502	6.3
SH0421	Phosphatases	299	2369	7.9
SH2821	T-Cell Activation	240	1265	5.3
SH0521	Tumor Suppressors	73	443	6.1
SH2121	Ubiquitin Ligases (E1, E2, E3)	203	1025	5.0

*The MISSION production and bio-informatics team constantly reviews and quality controls clones available for a gene family set. These numbers are very close to the actual number that will be shipped, but each researcher will receive a final plate map indicating the location and exact TRCN clone numbers.

Lentiviral Plasmid Vector pLKO.1-puro Features

Name	Description
U6	U6 Promoter
cppt	Central polypurine tract
hPGK	Human phosphoglycerate kinase eukaryotic promoter
puroR	Puromycin resistance gene for mammalian selection
SIN/3' LTR	3' self inactivating long terminal repeat
f1 ori	f1 origin of replication
ampR	Ampicillin resistance gene for bacterial selection
pUC ori	pUC origin of replication
5' LTR	5' long terminal repeat
Psi	RNA packaging signal
RRE	Rev response element



Control Selection Table

Sigma's recommended controls for any shRNA experiment are closely aligned with the controls suggested in the *Nature Cell Biology* editorial.⁴

Recommended Control	Objective
Negative Control: Untreated Cells	Untreated cells will provide a reference point for comparing all other samples.
Negative Control: Transfection with empty vector, containing no shRNA insert	MISSION pLKO.1-puro Control Vector, Catalog Number SHC001 The empty vector, pLKO.1-puro, is a useful negative control that will not activate the RNAi pathway because it does not contain an shRNA insert. It will allow for observation of cellular effects of the transfection process and the delivery of the lentiviral vector. Cells transfected with the empty vector provide a useful reference point for comparing specific knockdown.
Negative Control: Transfection with non-targeting shRNA	MISSION Non-Target shRNA Control Vector, Catalog Number SHC002 This non-targeting shRNA vector is a useful negative control that will activate RISC and the RNAi pathway, but does not target any human or mouse genes. The short-hairpin sequence contains 5 base pair mismatches to any known human or mouse gene. This allows for examination of the effects of shRNA transfection on gene expression. Cells transfected with the non-target shRNA vector will also provide a useful reference for interpretation of knockdown.
Positive Control: Transfection with positive reporter vector	MISSION TurboGFP Control Vector, Catalog Number SHC003 This vector is a useful positive control for measuring transfection efficiency and optimizing shRNA delivery. The TurboGFP Control vector consists of the lentiviral backbone vector, pLKO.1-puro, containing a gene encoding TurboGFP, driven by the CMV promoter. Transfection of this vector provides fast visual confirmation of successful transfection and delivery.
Positive Control: Transfection with shRNA targeting reporter vector	MISSION TurboGFP shRNA Control Vector, Catalog Number SHC004 The TurboGFP shRNA vector consists of the pLKO.1-Puro vector, containing shRNA that targets TurboGFP, and can be used as a positive control to quickly visualize knockdown. This TurboGFP shRNA Control Vector has been experimentally shown to reduce GFP expression by 99.6% in HEK 293T cells after 24 hours. Because this vector targets TurboGFP, and it does not target any human or mouse genes, it can also be used as a negative non-target control in shRNA experiments

Cell Type Table

The cell types listed below have been successfully infected by pLKO.1-puro based shRNA constructs

Cell lines, human	Cell Type	Cell lines, human	Cell Type	Primary cells human	Cell Type
HEK293	embryonic kidney cells	A431	epidermal carcinoma	dendritic	immature dendritic
HeLa	cervical adenocarcinoma	THP1	monocytic	T-cells	lymphocytes
A549	lung adenocarcinoma	RAW264.7	macrophage	epithelial	prostate
H1299	lung carcinoma	SH-SY5Y	brain neuroblastoma	fibroblasts	primary mammary
HT29-D4	colon carcinoma	HCN-1A	brain cortical neuron	Primary cells, other species	Cell Type
HepG2	hepatocellular carcinoma	SupT1	T-cells	ECS	mouse embryonic stem cells
HCT116	colon carcinoma	BJ-TERT	diploid fibroblasts	fibroblasts	mouse embryonic fibroblasts
MCF7	breast carcinoma	Cell lines, mouse	Cell Type	MC3T3-E1	mouse bone marrow derived
MCF10A	breast carcinoma	NIH3T3	fibroblast	molar mesenchymal	mouse embryonic mesenchymal
Panc-1	pancreatic epithelioid carcinoma	Primary cells, human	Cell Type	cardiomyocytes	rat neonatal cardiomyocytes
PC3	prostate carcinoma	astrocytes	normal		
DU145	prostate carcinoma	C3H10T1/2	mesenchymal		

References

1. Stewart, S.A., et al., Lentivirus-delivered stable gene silencing by RNAi in primary cells, *RNA*, **9**, 493-501 (2003).
2. Zufferey R, et al., Multiply attenuated lentiviral vector achieves efficient gene delivery *in vivo*, *Nat. Biotechnol.* **15**, 871-85 (1997).
3. Zufferey R, et al., Self-inactivating lentivirus vector for safe and efficient *in vivo* gene delivery, *J Virol.*, **72**, 9873-80 (1998).
4. Whither RNAi? *Nature Cell Biology*, **5**, 489-490 (2003).

Reviews Indicating the Importance of Each of the Gene Family Sets-

Apoptosis Pathway

1. Krysko, D.V., *et. al.* Apoptosis and necrosis: detection, discrimination and phagocytosis. *Methods*, **44**, 205-21 (2008).
2. Howley, B. and Fearnhead, H.O., Caspases as therapeutic targets. *J. Cell Mol. Med.*, Feb 24 [Epub ahead of print] (2008)
3. Logue, S.E. and Martin, S.J.. Caspase activation cascades in apoptosis. *Biochem. Soc. Trans.* **36 (Pt 1)**, 1-9 (2008).

B Cell Activation

1. Tolar, P., *et. al.* Viewing the antigen-induced initiation of B-cell activation in living cells. *Immunol Rev.*, **221**, 64-76 (2008).
2. Youinou, P., B cell conducts the lymphocyte orchestra. *J. Autoimmun.*, **28**, 143-51. (2007).

Cell Adhesion

1. Ebnet, K., Organization of multiprotein complexes at cell-cell junctions. *Histochem. Cell Biol.*, Mar 26 [Epub ahead of print] (2008).
2. Basson, M.D., An intracellular signal pathway that regulates cancer cell adhesion in response to extracellular forces. *Cancer Res.*, **68**, 2-4 (2008).
3. Mousa, S.A., Cell adhesion molecules: potential therapeutic & diagnostic implications. *Mol. Biotechnol.*, **38**, 33-40. (2008).

Cytokine and Chemokine Receptors

1. Callewaere, C, *et. al.* Chemokines and chemokine receptors in the brain: implication in neuroendocrine regulation. *J. Mol. Endocrinol.*, **38**, 355-63 (2007)
2. Allen, S.J., *et. al.* Chemokine: receptor structure, interactions, and antagonism. *Annu. Rev. Immunol.*; **25**, 787-820 (2007).
3. Zlotnik, A., *et. al.* The chemokine and chemokine receptor superfamilies and their molecular evolution. *Genome Biol.*; **7**, 243 (2006).
4. Mantovani, A., *et. al.* Regulatory pathways in inflammation. *Autoimmun. Rev.*, **7**, 8-11 (2007).

Cytokines and Chemokines

1. Anderson, P. Post-transcriptional control of cytokine production. *Nat. Immunol.*, **9**, 353-9 (2008).
2. Tayal, V. and Kalra, B.S., Cytokines and anti-cytokines as therapeutics--an update. *Eur. J. Pharmacol.*, **579**, 1-12 (2008).

Cytoskeleton

1. Dalby, M.J. and Yarwood, S.J., Analysis of focal adhesions and cytoskeleton by custom microarray. *Methods Mol. Biol.*, **370**, 121-34 (2007).
2. Dustin, M.L., Cell adhesion molecules and actin cytoskeleton at immune synapses and kinapses. *Curr. Opin. Cell Biol.*, **19**, 529-33 (2007).

DNA Repair Pathway

1. Hinkal, G. and Donehower, L.A., How does suppression of IGF-1 signaling by DNA damage affect aging and longevity? *Mech. Ageing Dev.*, **129**, 243-53 (2008).
2. Hakem, R., DNA-damage repair; the good, the bad, and the ugly. *EMBO J.*, **27**, 589-605 (2008).
3. Harper, J.W. and Elledge, S.J., The DNA damage response: ten years after. *Mol. Cell.*, **28**, 739-45 (2007).

DUBS - Ubiquitin Hydrolyases

1. Nicholson, B, *et. al.* Deubiquitinating enzymes as novel anticancer targets. *Future Oncol.*, **3**, 191-9 (2007).
2. Millard, S.M. and Wood, S.A., Riding the DUBway: regulation of protein trafficking by deubiquitylating enzymes. *J. Cell Biol.*, **173**, 463-8 (2006).
3. Amerik, A.Y. and Hochstrasser. M., Mechanism and function of deubiquitinating enzymes. *Biochim. Biophys. Acta*, **1695**, 189-207 (2004).

Epigenetic Regulators

1. Esteller, M., Epigenetics in cancer. *N. Engl. J. Med.*, **358**, 1148-59. Review (2008).
2. Grønbaek, K., *et. al.* Epigenetic changes in cancer. *APMIS*, **115**, 1039-59 (2007).

Extracellular Matrix

1. Rees, M.D., *et. al.* Oxidative damage to extracellular matrix and its role in human pathologies. *Free Radic. Biol. Med.*, Apr 8 (2008). [Epub ahead of print]
2. Adair-Kirk, T.L. and Senior, R.M., Fragments of extracellular matrix as mediators of inflammation. *Int. J. Biochem. Cell Biol.*, **40**, 1101-10 (2008).
3. Daley, W.P., *et. al.* Extracellular matrix dynamics in development and regenerative medicine. *J. Cell Sci.*, **121(Pt 3)**, 255-64 (2008).

G-Protein-Coupled Receptors:

1. Thompson, M.D., *et. al.* G protein-coupled receptors disrupted in human genetic disease. *Methods Mol. Biol.*; **448**, 109-37 (2008).
2. Milligan, G., New aspects of G-protein-coupled receptor signalling and regulation. *Trends Endocrinol. Metab.*, **9**, 13-9 (1998).

Helicases

1. Ha, T., Need for speed: mechanical regulation of a replicative helicase. *Cell*, **129**, 1249-50 (2007).
2. Singleton, M.R., et al., Structure and mechanism of helicases and nucleic acid translocases. *Annu. Rev. Biochem.*, **76**, 23-50 (2007).
3. Xi, X.G., Helicases as antiviral and anticancer drug targets. *Curr. Med. Chem.*, **14**, 883-915 (2007).

Ion Channels

1. Cannon, S.C., Physiologic principles underlying ion channelopathies. *Neurotherapeutics*, **4**, 174-83 (2007).

JAK-STAT Pathway

1. Murray, P.J., The JAK-STAT signaling pathway: input and output integration. *J. Immunol.*, **178**, 2623-9 (2007).
2. O'Sullivan, L.A., et al. Cytokine receptor signaling through the Jak-Stat-Socs pathway in disease. *Mol. Immunol.*, **44**, 2497-506 (2007).

Kinases

1. Gomase, V.S., et al., *Curr. Drug Metab.*, **9**, 255-8 (2008).

Nuclear Hormone Receptors

1. Kininis, M. and Kraus, W.L., A global view of transcriptional regulation by nuclear receptors: gene expression, factor localization, and DNA sequence analysis. *Nucl. Recept. Signal*, **6**, e005 (2008).

p53 Pathway

1. Bose, I. And Ghosh, B., The p53-MDM2 network: from oscillations to apoptosis. *J. Biosci.*, **32**, 991-7 (2007).
2. Efeyan, A. and Serrano, M., p53: guardian of the genome and policeman of the oncogenes. *Cell Cycle*, **6**, 1006-10 (2007).
3. Kastan, M.B., Wild-type p53: tumors can't stand it. *Cell*, **128**, 837-40 (2007).

Phosphatases

1. Hendriks, W.J., et al. Protein tyrosine phosphatases: functional inferences from mouse models and human diseases. *FEBS J.*, **275**, 816-30 (2008).
2. Tremblay, M.L. and Giguère, V., Phosphatases at the heart of FoxO metabolic control. *1: Cell Metab.*, **7**, 101-3 (2008).
3. Heideker, J., et al. Phosphatases, DNA damage checkpoints and checkpoint deactivation. *Cell Cycle*, **6**, 3058-64 (2007).
4. Sawyer, T.K., et al. Protein phosphorylation and signal transduction modulation: chemistry perspectives for small-molecule drug discovery. *Med. Chem.*, **1**, 293-319 (2005).

T Cell Activation

1. Won, J. and Lee, GH., T-cell-targeted signaling inhibitors. *Int. Rev. Immunol.*, **27**, 19-41 (2008).
2. Brenner, D., et al. Concepts of activated T cell death. *Crit. Rev. Oncol. Hematol.*, **66**, 52-64 (2008).
3. Seminario, M.C. and Bunnell, S.C., Signal initiation in T-cell receptor microclusters. *Immunol. Rev.*, **221**, 90-106 (2008).
4. Lämmermann, T, and Sixt, M. The microanatomy of T-cell responses. *Immunol. Rev.*, **221**, 26-43 (2008).

Tumor Suppressors

1. Vatteemi, E. and Claudio, P.P., Tumor suppressor genes as cancer therapeutics. *Drug News Perspect*, **20**, 511-20 (2007).
2. Berger, J.C. et al. Metastasis suppressor genes: from gene identification to protein function and regulation. *Cancer Biol. Ther.*, **4**, 805-12 (2005).

Ubiquitin Ligases (E1, E2, E3)

1. Cardozo, T. and Pagano, M., Wrenches in the works: drug discovery targeting the SCF ubiquitin ligase and APC/C complexes. *BMC Biochem.*, **8 Suppl 1**, S9 (2007).
2. Newton, K. and Vucic, D., Ubiquitin ligases in cancer: ushers for degradation. *Cancer Invest.*, **25**, 502-13 (2007).
3. Sun, Y. Overview of approaches for screening for ubiquitin ligase inhibitors. *Methods Enzymol.*, **399**, 654-63 (2005).
4. Hershko, A. The ubiquitin system for protein degradation and some of its roles in the control of the cell division cycle. *Cell Death Differ.*, **12**, 1191-7 (2005).

MISSION is a registered trademark of Sigma-Aldrich Biotechnology LP and Sigma-Aldrich Co.
TurboGFP is a trademark of Evrogen Co.

Limited Use Licenses

Sigma has acquired necessary key licenses for lentiviral systems and RNAi and provides freedom to operate under our label license for relevant purchased products. Because Sigma actively evaluates this rapidly evolving intellectual property space, please visit www.sigma.com/shrna for up-to-date information on current licenses for the MISSION® shRNA collections.

Use of this product for Commercial Purposes requires a license from Sigma-Aldrich Corporation. The purchase of this product conveys to the buyer the nontransferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party, or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. Commercial Purposes means any activity by a party for consideration, but excludes not-for-profit core facilities providing services within their own research institutions at cost. Core facilities are invited to join Sigma-Aldrich's RNAi Partnership Program. Details of Sigma-Aldrich's RNAi Partnership Program can be found at www.sigma.com/rpp.

The MISSION shRNA Library of The RNAi Consortium is produced and distributed under license from the Massachusetts Institute of Technology.

Licensed under Carnegie Institution US Patent 6,506,559 and Massachusetts Institute of Technology and for laboratory and commercial research use only.

This product is licensed under U.S. Pat. Nos. 5,817,491; 5,591,624; 5,716,832; 6,312,682; 6,669,936; 6,235,522; 6,924,123 and foreign equivalents from Oxford BioMedica (UK) Ltd., Oxford, UK, and is provided for use in academic and commercial *in vitro* and *in vivo* research for elucidating gene function, and for validating potential gene products and pathways for drug discovery and development, but excludes any use of LentiVector® technology for: creating transgenic birds for the purpose of producing useful or valuable proteins in the eggs of such transgenic birds, the delivery of gene therapies, and for commercial production of therapeutic, diagnostic or other commercial products not intended for research use where such products do not consist of or incorporate a lentiviral vector. Information about licenses for commercial uses excluded under this license is available from Oxford BioMedica (UK), Ltd., Medawar Center, Oxford Science Park, Oxford OX4 4GA UK enquiries@oxfordbiomedica.co.uk or BioMedica Inc., 11622 El Camino Real #100, San Diego CA 92130-2049 USA. LentiVector is a registered US and European Community trademark of Oxford BioMedica plc.

This product is licensed under agreement with Benitec Australia Ltd. and CSIRO as co-owners of U.S. Pat. No. 6,573,099 and foreign counterparts, for use in research to understand, diagnose, monitor, treat and prevent human diseases and disorders, including the use of animals for such research use, except that use of ddRNAi as a therapeutic agent or as a method of disease treatment, prevention, diagnosis or for disease monitoring is excluded. Information regarding licenses to these patents for use of ddRNAi as a therapeutic agent or for other uses excluded under this license is available from Benitec at licensing@benitec.com. Information about licenses for the use of ddRNAi in other fields, is available from CSIRO at pi.csiro.au/RNAi.

This product is licensed under European Patents 1144623, 1214945 and foreign equivalents from Alnylam Pharmaceuticals, Inc., Cambridge, USA and is provided only for use in academic and commercial research whose purpose is to elucidate gene function, including research to validate potential gene products and pathways for drug discovery and development and to screen non-siRNA based compounds (but excluding the evaluation or characterization of this product as the potential basis for a siRNA-based drug) and not for any other commercial purposes. Information about licenses for commercial use (including discovery and development of siRNA-based drugs) is available from Alnylam Pharmaceuticals, Inc., 300 Third Street, Cambridge, MA 02142, USA.

This product (based upon the lentikat system) is sub-licensed from Invitrogen Corporation under U.S. Patent Nos. 5,686,279, 5,834,256, 5,858,740; 5,994,136; 6,013,516; 6,051,427, 6,165,782, and 6,218,187 and corresponding patents and applications in other countries for internal research purposes only. Use of this technology for gene therapy applications or bioprocessing other than for nonhuman research use requires a license from Cell Genesys, Inc. Please contact Cell Genesys, Inc. at 342 Lakeside Drive, Foster City, California 94404. Use of this technology to make or sell products or offer services for consideration in the research market requires a license from Invitrogen Corporation, 1600 Faraday Ave., Carlsbad, CA 92008.

This product is for non-clinical research use only. It is not to be used for commercial purposes. Use of this product to produce products for sale or for diagnostic, therapeutic or high throughput drug discovery purposes (the screening of more than 10,000 compounds per day) is prohibited. This product is sold under license from Invitrogen Corporation. In order to obtain a license to use this product for these commercial purposes, contact The Regents of the University of California. This product or the use of this product is covered by U.S. Patent No. 5,624,803 owned by The Regents of the University of California.

BR,CB,PHC 10/09-1

Sigma brand products are sold through Sigma-Aldrich, Inc.
Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications.
Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply.
Please see reverse side of the invoice or packing slip.