

SUPERKILLERTRAIL®

Storage and Dilution Buffer

Storage and dilution buffer for use with the following product:
SUPERKILLERTRAIL®, Soluble (human) (recombinant) (Prod. No. ALX-201-115)

Useful for prediluting the above mentioned product before preparing aliquots or performing experiments. Predilution with SUPERKILLERTRAIL® Storage and Dilution Buffer allows for easier handling of small working concentrations (sometimes only 0.1µl of the SUPERKILLERTRAIL® stock solution is necessary because of their high activity). Predilution with SUPERKILLERTRAIL® Storage and Dilution Buffer instead of medium or PBS prevents loss of biological activity. Can also be used as negative control.

Ordering Information

[Order Online »](#)

ENZ-PRT317-0500	500µl
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.

Long Term Storage -80°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Formulation Liquid. 20 mM Hepes, 120 mM NaCl, 0.01% Tween20, 1% Sucrose, 1 mM DTT, pH = 7.9.

Purity Detail Sterile filtered (0.22 µm).

Technical Info / Product Notes Prepare a 1:5 dilution of SUPERKILLERTRAIL[®] with SUPERKILLERTRAIL[®] Storage and Dilution Buffer. For immediate use, keep prediluted SUPERKILLERTRAIL[®] on ice until adding to cells.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)