

# KillerTRAIL™ Storage and Dilution Buffer

Storage and dilution buffer for use with the following product:

*KillerTRAIL™* Protein (soluble) (human), (recombinant) (Prod. No. ALX-201-073)

Useful for prediluting the above mentioned product before preparing aliquots or performing experiments. Predilution with *KillerTRAIL™* Storage and Dilution Buffer allows for easier handling of small working concentrations (sometimes only 0.1µl of the *KillerTRAIL™* stock solution is necessary because of their high activity. Predilution with *KillerTRAIL™* 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-PRT318-0500	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, 300 mM NaCl, 0.01% Tween20, 1% Sucrose, 1 mM DTT, pH = 7.4.
Purity Detail	Sterile filtered (0.22 µm).
Technical Info / Product Notes	Prepare a 1:5 dilution of <i>KillerTRAIL</i> <sup>™</sup> with <i>KillerTRAIL</i> <sup>™</sup> Storage and Dilution Buffer. For immediate use, keep prediluted <i>KillerTRAIL</i> <sup>™</sup> 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)