## KillerTRAIL™ Storage and Dilution Buffer

Storage and dilution buffer for use with the following product: *Killer*TRAIL™ Protein (soluble) (human), (recombinant) (Prod. No. ALX-201-073)

Useful for prediluting the above mentioned product before preparing aliquots or performing experiments. Predilution with *Killer*TRAIL™ 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 *Killer*TRAIL™ 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 KillerTRAIL™ with KillerTRAIL™

> Storage and Dilution Buffer. For immediate use, keep prediluted KillerTRAIL™ on ice until adding to cells.

Last modified: May 29, 2024



info-

eu@enzolifesciences.com