SUPERKILLERTRAIL® Storage and Dilution Buffer

Storage and dilution buffer for use with the following product: ${\sf SUPERKILLERTRAIL}^{\circledR} \ , \ {\sf Soluble} \ (human) \ (recombinant) \ (Prod. \ No. \ ALX-201-115)$

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).

Prepare a 1:5 dilution of SUPERKILLERTRAIL® with **Technical Info / Product Notes**

> SUPERKILLERTRAIL® Storage and Dilution Buffer. For immediate use, keep prediluted SUPERKILLERTRAIL® on

ice until adding to cells.

Last modified: May 29, 2024



info-

eu@enzolifesciences.com