

# Bax (human) (58-75), (cell permeable)

For increased cell permeability the peptide was covalently linked to the 10 aa recognized by the TAT transporter.

**Manuals, SDS & CofA**

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended.
<b>Handling</b>	Keep on ice at all times. Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
<b>Short Term Storage</b>	-20°C
<b>Long Term Storage</b>	-80°C
<b>Shipping</b>	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	BCL2L4 (58-75)
<b>Concentration</b>	10mM
<b>Formula</b>	$C_{144}H_{260}N_{58}O_{44}S_2$
<b>Formulation</b>	Liquid. In sterile PBS.
<b>MW</b>	3572.2
<b>Purity</b>	≥95% (HPLC)
<b>Quantity</b>	50µl
<b>Sequence</b>	GRKKRRQRRR-GG-KLSECLKRIGDELDSNME-amide

Last modified: May 29, 2024



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)