

# DT-3

## Peptide-based inhibitor of protein kinase G

Membrane-permeant peptide-based inhibitor of cyclic GMP-dependent protein kinase  $\alpha$  and  $\beta$  with nanomolar  $K_i$  values. It displays high selectivity for cyclic GMP-dependent protein kinase  $\alpha$  compared to cyclic AMP-dependent protein kinase with a ratio of approximately 20,000 fold (Dostmann *et al.* 2000).

*BLG-D052-025 (5 x 0.5 mg pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

### Ordering Information

[Order Online »](#)

BLG-D052-005	0.5mg
BLG-D052-025	5x0.5mg

### Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** PKG Inhibitor Peptide DT-3 . trifluoroacetate salt

**Appearance** Crystallized or lyophilized trifluoroacetate salt.

**Formula**  $C_{152}H_{259}N_{53}O_{27}S \cdot (C_2F_3O_2)_{16}$

**MW** 3293.1 (free acid)

**Purity** > 95% HPLC

**Quantity** 0.5mg~0.16µmol

**Sequence** RQIKIWFQNRRMKWKK-LRKKKKKKH-amide

**Solubility** Soluble in water (at least 10mM) or DMSO.

**Technical Info / Product Notes** For the Original Manufacturer's data sheet please [click here](#).



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)