

# W45

## Peptide-based, membrane-impermeant inhibitor of protein kinase G I $\alpha$ and I $\beta$

W45 is a potent peptide-based, membrane-impermeant inhibitor of cyclic GMP-dependent protein kinase I $\alpha$  and I $\beta$ . It displays selectivity for cyclic GMP-dependent protein kinase I $\alpha$  compared to cyclic AMP-dependent protein kinase with a ratio of approximately 600 fold (Dostmann *et al.* 2000).

Fusion of W45 to membrane translocation sequences from HIV-1 Tat protein or *Drosophila Antennapedia* homeodomain results in the membrane-permeant fusion peptides DT-2 (Prod. No. BLG-D051) and DT-3 (Prod. No. BLG-D052), respectively. Since the highly charged W45 cannot cross cell membranes it can be used as a control in applications of DT-2, DT-3 and the proteolytically stable D-amino acid congener of DT-2, (D)-DT-2 (Prod. No. BLG-D110).

*BLG-W001-025 (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-W001-005	0.5mg
BLG-W001-025	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 W45 . trifluoroacetate salt

**Formula** C<sub>48</sub>H<sub>93</sub>N<sub>19</sub>O<sub>8</sub> (free peptide)

**MW** 1064.4 (free peptide)

**Purity** > 95% HPLC

**Reconstitution** Addition of 470µl water or buffer to a 0.5mg vial yields a 1mM solution.

**Sequence** L-R-K-K-K-K-K-H-amide (W45)

**Solubility** Soluble in water.

**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)