

(D)-DT-2

Potent peptide-based inhibitor of protein kinase G I ? and I ?

(D)-DT-2 is a peptide-based, membrane-permeant and proteolytically stable inhibitor of cyclic GMP-dependent protein kinase (PKG) I α and I β with nanomolar *in vitro* IC₅₀ values (Nickl et al. 2010, Gambaryan et al. 2011). It displays high selectivity for cyclic GMP-dependent protein kinase I α compared to cyclic AMP-dependent protein kinase with a ratio of approximately 15,000 fold (Nickl et al. 2010). (D)-DT-2 is the newly developed D-amino acid analogue of the peptidic PKG inhibitor DT-2 (Prod. No. BLG-D051). So far, the inhibitory potential of (D)-DT-2 in *in vivo* applications has not been fully characterized, but seems to be variable depending on the type of biosystem used. It is recommended to control the physiological effects of (D)-DT-2 in *in vivo* applications by PKG specific substrate phosphorylation.

BLG-D110-025 (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-D110-005	0.5mg
BLG-D110-025	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 (D)-DT-2 . trifluoroacetate salt

Formula $C_{122}H_{223}N_{53}O_{23}$ (free peptide)

MW 2800.5 (free peptide)

Purity > 95% HPLC

Sequence YGRKKRRQRRRPPLRKKKKKH-NH₂ (all-d)

Solubility Soluble in water (≥6mM).

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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com