

Boc-DON-Gln-Ile-Val-OMe

Selective and irreversible inhibitor of tissue transglutaminase [TG2; protein-glutamine γ -glutamyltransferase]. The reactive group of this inhibitor is a modified glutamic acid (6-diazo-5-oxonorleucin) [1;2].

Ordering Information

[Order Online »](#)

NZY-B003-010	10mg
--------------	------

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 Boc-L-(6-diazo-5-oxonorleucinyI)-Gln-Ile-Val-OMe

Appearance Yellow solid.

Formula $C_{28}H_{47}N_7O_9$

MW 625.7

Purity ≥95% (HPLC).

Solubility Predissolve in DMSO. Soluble (1.5mM) in 2% DMSO/aqueous buffer.

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