

L-TAT control peptide (FITC conjugate)

Control peptide for L-JNKI

Control peptide for L-JNKI. For more details see L-JNKI1 (Prod.No. BML-EI354).

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-168-009F-R050

FITC - 50µl

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	JNK Inhibitor 1 (L-stereoisomer) control peptide (FITC conjugate)
Concentration	1mM
Formula	$C_{80}H_{128}N_{34}O_{17}S$
Formulation	Liquid. In sterile PBS.
MW	1870.2
Purity	≥90% (HPLC)
Sequence	FITC-β-A-RKKRRQRRRA-amide



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-fr@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