

RK-682

VHR inhibitor

RK-682 is a novel tyrosine phosphatase inhibitor isolated from *Streptomyces sp.* It inhibits the dual specificity VHR phosphatase (IC_{50} =2.0 μ M) selectively over CD45 (IC_{50} =54 μ M), Cdc25B and PTP1B (IC_{50} ~42 μ M). RK-682 inhibits cell cycle progression of Ball-1 cells at the G₁/S transition and enhances ATP-induced long-term potentiation in guinea pig hippocampal CA1 neurons. Shows potent activity against HIV-1 protease.

Citations: 6

[View Online »](#)

Ordering Information

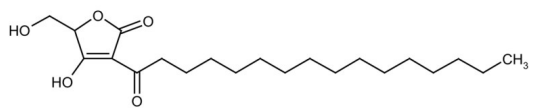
[Order Online »](#)

BML-PR112-0200

200 μ g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance	Pale yellow solid.
CAS	154639-24-4
Couple Target	Serine/threonine-protein phosphatase
Couple Type	Inhibitor
Formula	$C_{21}H_{36}O_5$
MW	368.5
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (25mg/ml), 100% ethanol (10mg/ml), methanol or dimethyl formamide.

Last modified: May 29, 2024



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