

# RK-682

## VHR inhibitor

RK-682 is a novel tyrosine phosphatase inhibitor isolated from *Streptomyces sp.* It inhibits the dual specificity VHR phosphatase ( $IC_{50}=2.0\mu M$ ) selectively over CD45 ( $IC_{50}=54\mu M$ ), Cdc25B and PTP1B ( $IC_{50}\sim 42\mu M$ ). RK-682 inhibits cell cycle progression of Ball-1 cells at the G<sub>1</sub>/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 »](#)

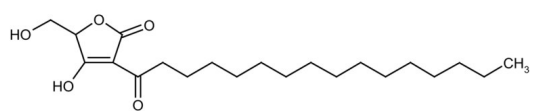
## Ordering Information

[Order Online »](#)

BML-PR112-1000	1mg
BML-PR112-0200	200µg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	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.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

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



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)