

RWJ-60475 (AM)₃

CD45 inhibitor

RWJ-60475-(AM)₃ is a cell-permeable acetoxymethyl (AM) ester of RWJ-60475. The AM esters confer cell membrane permeability and are hydrolyzed by intracellular esterases, trapping the active inhibitor within the cell. The IC₅₀ for the hydrolyzed, active inhibitor is 2 μM.

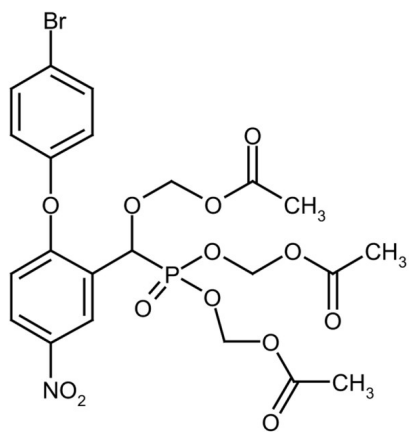
Ordering Information

[Order Online »](#)

BML-PR110-0001	1mg
----------------	-----

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. Make solutions fresh daily.
Handling	Protect from moisture.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance	Pale yellow-brown oil.
Couple Target	CD45, PTP
Couple Type	Inhibitor
Formula	$C_{22}H_{23}BrNO_{13}P$
MW	620.4
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO (25 mg/ml), acetone, or dimethylchloride.

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