

BMOV

Potent insulin mimetic

Organic vanadium complex which acts as a potent insulin mimic. It normalizes glucose and lipid values and restores food and fluid intake to control levels in diabetic rats without increasing insulin levels (0.37 mmol/kg). It attenuates hyperinsulinemia and hypertension in spontaneously hypertensive rats. Its mechanism of action includes inhibition of PTP1B as well as other tyrosine phosphatases. The drug is nontoxic *in vitro* at useful concentrations and displays no signs of toxicity in mice at high oral doses.

Citations: 2

[View Online »](#)

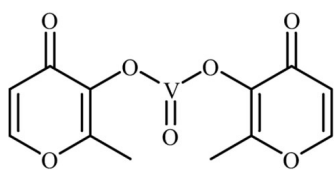
Ordering Information

[Order Online »](#)

BML-PR131-0050	50mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at room temperature. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Bis(maltolato)oxovanadium (IV)
Appearance	Black solid.
CAS	38213-69-3
Couple Target	PTP
Couple Type	Inhibitor
Formula	$C_{12}H_{10}O_7V$
MW	317.2
Purity	≥96% (TLC)
Solubility	Soluble in water (10mg/ml) and DMSO (15mg/ml).



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)