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

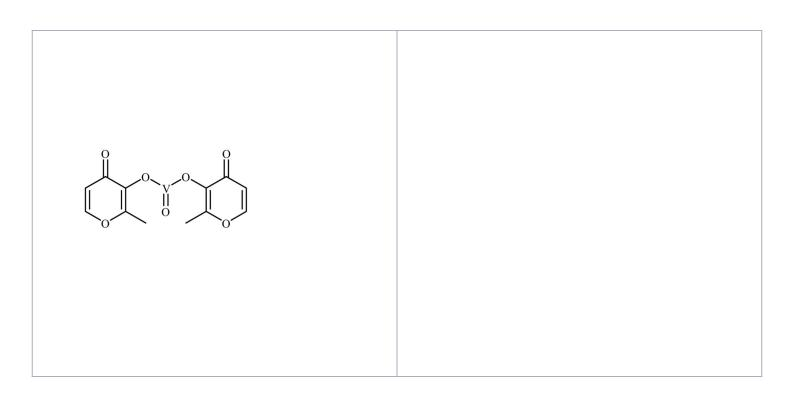
Order Online »

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)

Soluble in water (10mg/ml) and DMSO (15mg/ml).

Last modified: May 29, 2024