

Cantharidic acid

PP2A inhibitor

Potent and selective inhibitor of PP2A, $IC_{50}=53nM$ (IC_{50} for PP1=562nM). Useful for inhibiting PP1 and PP2A during purification of phosphorylated proteins.

Citations: 3

[View Online »](#)

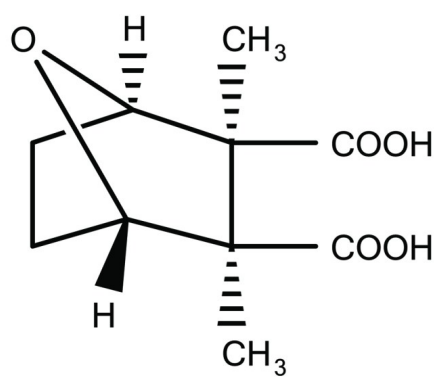
Ordering Information

[Order Online »](#)

| | |
|----------------|------|
| BML-PR106-0010 | 10mg |
| BML-PR106-0050 | 50mg |

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

| | |
|-------------------|--|
| Use/Stability | As indicated on product label or CoA when stored as recommended. Store, as supplied, at room temperature for up to 5 years. Store solutions at -20°C for up to 3 months. |
| Long Term Storage | Ambient |
| Shipping | Ambient Temperature |

Regulatory Status RUO - Research Use Only

Product Details

| | |
|---------------|--|
| Appearance | White solid. |
| CAS | 28874-45-5 |
| Couple Target | Serine/threonine-protein phosphatase |
| Couple Type | Inhibitor |
| Formula | $C_{10}H_{14}O_5$ |
| MW | 214.2 |
| MeltingPoint | 216-218°C |
| Purity | ≥98% (TLC) |
| Solubility | Soluble in DMSO (25mg/ml) or ethanol (10mg/ml warm). |



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-
fr@enzolifesciences.com](mailto:info-fr@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)