

Tranylcypromine

Demethylase inhibitor

Inhibits prostacyclin synthase. Monoamine oxidase inhibitor. Effective small molecule inhibitor of histone demethylation. Inhibits the demethylase BHC110/LSD1 ($IC_{50} < 2\mu M$). When combined with a GSK-3 inhibitor, it causes reprogramming of human primary keratinocyte transduced with Oct4 and Klf4.

Citations: 8

[View Online »](#)

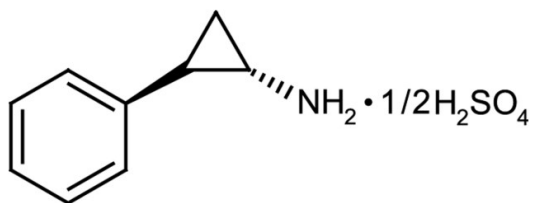
Ordering Information

[Order Online »](#)

BML-EI217-0001	1g
BML-EI217-0005	5g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature – Dangerous Good

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Trans(±)-2-phenylcyclopropanamine, SKF trans-355, Tranlylcypromine hemisulfate

Appearance White solid.

CAS 13492-01-8

Couple Target Demethylase

Couple Type Inhibitor

Formula $C_9H_{11}N \cdot 1/2H_2SO_4$

MW 182.2

Purity ≥98% (HPLC)

Solubility Soluble in water (25mg/ml).

Last modified: May 29, 2024



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)