

(S)-Mephenytoin

Cytochrome P450 substrate

(S)-Mephenytoin is a substrate of CYP2C19 for the analysis of cytochrome P450 metabolism and polymorphism. (S)-Mephenytoin is an anti-convulsant and anti-epileptic.

Citations: 1

[View Online »](#)

Ordering Information

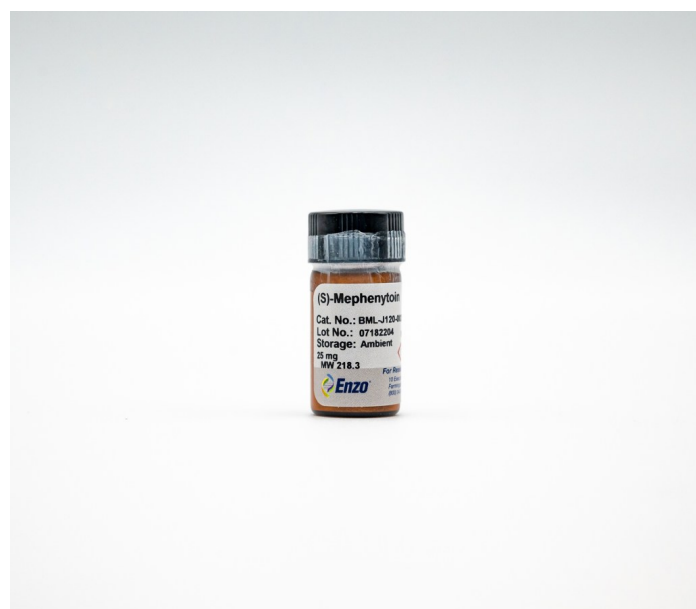
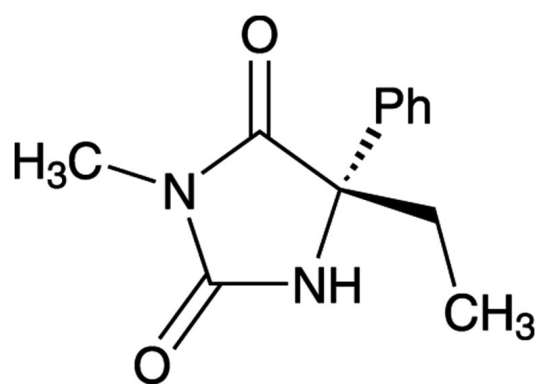
[Order Online »](#)

BML-J120-0005	5mg
BML-J120-0025	25mg

Manuals, SDS & CofA

[View Online »](#)

- Highly pure compound used as a substrate for the screening of genetic defects in CYP2C19



Handling & Storage

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

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name (+)-Mephencytoin, (S)-5-ethyl-3-methyl-5-phenylhydantoin

Appearance White solid.

CAS 70989-04-7

Couple Target Cytochrome P450

Couple Type Substrate

Formula $C_{12}H_{14}N_2O_2$

Identity Identity determined by 1H-NMR.

MW 218.3

Purity ≥98% (HPLC)

Solubility Soluble in DMSO or 100% ethanol.

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)