

Ethacrynic acid

GST inhibitor

Inhibits LTC₄ biosynthesis by inhibition of glutathione S-transferase (GST).
Inhibits 15-hydroxyprostaglandin dehydrogenase.

Citations: 3

[View Online »](#)

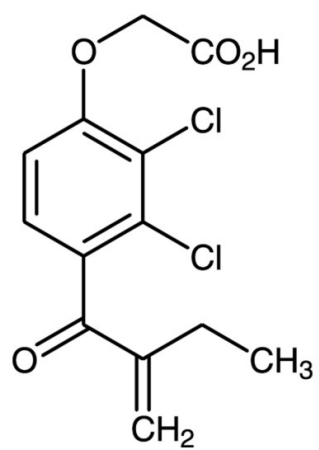
Ordering Information

[Order Online »](#)

BML-EI128-0001	1g
----------------	----

Manuals, SDS & CofA

[View Online »](#)



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 2,3-Dichloro-4-(2-methylene-1-oxobutyl)phenoxyacetic acid

Appearance Off-white to white.

CAS 58-54-8

Couple Target GST

Couple Type Inhibitor

MW 303.1

Purity ≥98%

Solubility Soluble in 100% ethanol (35mg/ml) or DMSO (35mg/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)