

α -Hydroxytriazolam

An active but short-lived metabolite of triazolam

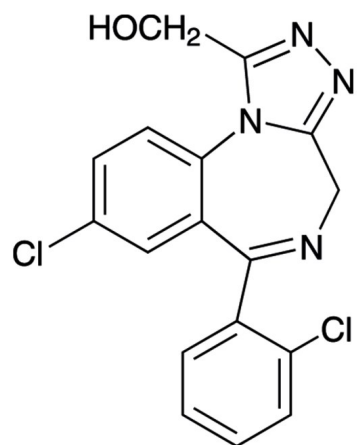
Metabolite of triazolam.

Citations: 4

[View Online »](#)

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

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 8-Chloro-6-(o-chlorophenyl)-4H-s-Triazolo[4,3-a][1,4]-benzodiazepine-1-methanol

Appearance White solid.

CAS 37115-45-0

Formula $C_{17}H_{12}Cl_2N_4O$

Identity Determined by 1H-NMR and MS.

MW 359.2

Purity ≥96% (TLC)

Solubility Soluble in DMSO (5mg/ml warm) or methanol (1mg/ml warm).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-de@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com