

Zolpidem Metabolite I

Major metabolite of zolpidem

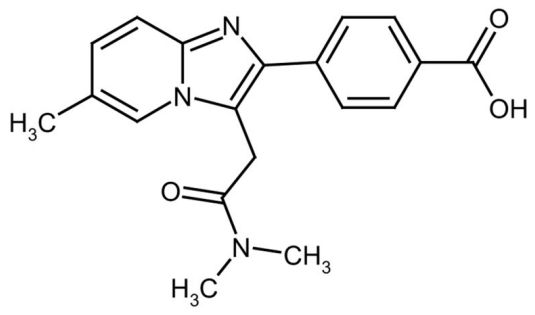
Zolpidem Metabolite I is the major metabolite of zolpidem. Zolpidem is a non-benzodiazepine compound that acts as an agonist to the GABA-A γ 1 subunit, which has been shown to treat insomnia. Intake of zolpidem can be detected by quantification of the metabolites produced. Zolpidem metabolite I shows two absorption peaks in methanol. It emits blue fluorescence upon exciting at 303 nm in water ($\lambda_{em} = 373$ nm).

Citations: 1

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Hygroscopic. Protect from moisture. Keep cool and dry.
Long Term Storage	-20°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	4-[3-[2-(Dimethylamino)-2-oxoethyl]-6-methylimidazo[1,2-a]pyridin-2-yl] benzoic acid, Zolpidem carboxylic acid
Appearance	Off-white solid.
CAS	109461-65-6
Couple Target	GABA receptor
Excitation Maximum	247nm and 303nm
Formula	$C_{19}H_{19}N_3O_3$
Identity	Determined by MS and 1H-NMR.
MW	337.37
Purity	≥98% (HPLC)
Solubility	Soluble in methanol, DMF and DMSO. Slightly soluble in water and ethyl acetate.

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-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@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