

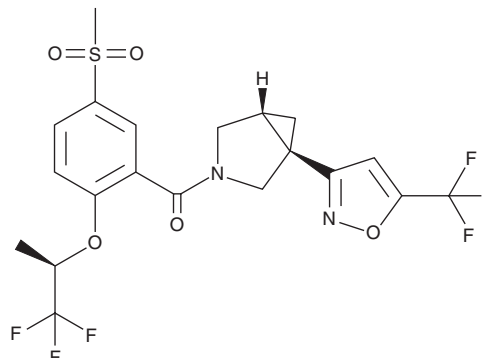
PRODUCT INFORMATION



Iclepertin

Item No. 41729

CAS Registry No.: 1421936-85-7
Formal Name: [5-(methylsulfonyl)-2-[(1R)-2,2,2-trifluoro-1-methylethoxy]phenyl][(1R,5R)-1-[5-(trifluoromethyl)-3-isoxazolyl]-3-azabicyclo[3.1.0]hex-3-yl]-methanone
Synonym: BI-425809
MF: C₂₀H₁₈F₆N₂O₅S
FW: 512.4
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Iclepertin is supplied as a solid. A stock solution may be made by dissolving the iclepertin in the solvent of choice, which should be purged with an inert gas. Iclepertin is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

Iclepertin is an inhibitor of glycine transporter 1 (GlyT1; IC₅₀s = 5 and 5.2 nM for the human and rat transporter, respectively).¹ It is selective for GlyT1 over GlyT2 (IC₅₀ = >10 μM for the human transporter). Iclepertin (0.2, 0.6, or 2 mg/kg) increases cerebrospinal fluid (CSF) levels of glycine in rats. It improves learning and memory in the social recognition test in rats when administered at doses of 0.2, 0.6, or 1.8 mg/kg.² Iclepertin (0.5, 1.5, or 4.5 mg/kg) reverses MK-801-induced working memory deficits in the T-maze test in mice.

References

1. Rosenbrock, H., Desch, M., Kleiner, O., *et al.* Evaluation of pharmacokinetics and pharmacodynamics of BI 425809, a novel GlyT1 inhibitor: Translational studies. *Clin. Transl. Sci.* **11**(6), 616-623 (2018).
2. Rosenbrock, H., Dorner-Ciossek, C., Giovannini, R., *et al.* Effects of the glycine transporter-1 inhibitor iclepertin (BI 425809) on sensory processing, neural network function, and cognition in animal models related to schizophrenia. *J. Pharmacol. Exp. Ther.* **382**(2), 223-232 (2022).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 01/27/2025

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM