

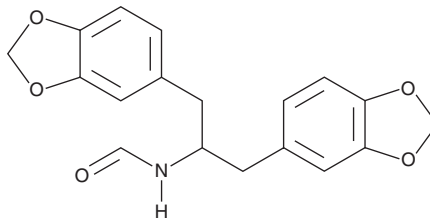
PRODUCT INFORMATION



ECSI#6

Item No. 42250

CAS Registry No.: 860456-56-0
Formal Name: N-[2-(1,3-benzodioxol-5-yl)-1-(1,3-benzodioxol-5-ylmethyl)ethyl]-formamide
MF: $C_{18}H_{17}NO_5$
FW: 327.3
Purity: $\geq 95\%$
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

ECSI#6 is supplied as a solid. A stock solution may be made by dissolving the ECSI#6 in the solvent of choice, which should be purged with an inert gas. ECSI#6 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile and water.

Description

ECSI#6 is an inhibitor of the serotonin transporter (SERT), dopamine transporter (DAT), and norepinephrine transporter (NET).¹ It inhibits serotonin, dopamine, and norepinephrine uptake in HEK293 cells expressing human SERT, DAT, and NET (IC_{50} s = 8, 48.2, and 17 μM , respectively).

Reference

1. Pifl, C., Nagy, G., Berényi, S., *et al.* Pharmacological characterization of ecstasy synthesis byproducts with recombinant human monoamine transporters. *J. Pharmacol. Exp. Ther.* **14**(1), 346-354 (2005).

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, 10/25/2024

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