

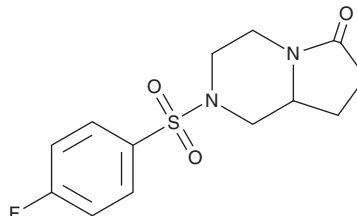
PRODUCT INFORMATION



Unifiram

Item No. 34209

CAS Registry No.: 272786-64-8
Formal Name: 2-[(4-fluorophenyl)sulfonyl]hexahydro-
pyrrolo[1,2-a]pyrazin-6(2H)-one
Synonym: DM232
MF: C₁₃H₁₅FN₂O₃S
FW: 298.3
Purity: ≥98%
UV/Vis.: λ_{max}: 223 nm
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

Unifiram is supplied as a solid. A stock solution may be made by dissolving the unifiram in the solvent of choice, which should be purged with an inert gas. Unifiram is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of unifiram in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Unifiram is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, unifiram should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Unifiram has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Unifiram is a nootropic agent.¹ It increases acetylcholine (ACh) release in the rat cerebral cortex *in vivo* and induces a long-lasting increase in the amplitude of field excitatory postsynaptic potentials (fEPSPs) in the rat hippocampal CA1 region (EC₅₀ = 27 nM). It does not bind to serotonin (5-HT), dopamine, muscarinic, nicotinic, adrenergic, glutamate, histamine, opioid, or GABA receptors at 1 μM. Unifiram (0.1 mg/kg) improves memory in non-memory-impaired rats in a social learning test.² It also prevents memory deficits induced by the anticholinergic agent scopolamine, nicotinic receptor antagonist mecamylamine, GABA_B receptor agonist baclofen, or α₂-adrenergic receptor agonist clonidine in the passive avoidance test in mice when administered at a dose of 0.01 mg/kg and prevents memory deficits induced by the AMPA/kainate glutamate receptor antagonist NBQX at 0.1 mg/kg.¹

References

1. Romanelli, M.N., Galeotti, N., Ghelardini, C., *et al.* Pharmacological characterization of DM232 (unifiram) and DM235 (sunifiram), new potent cognition enhancers. *CNS Drug Rev.* **12**(1), 39-52 (2006).
2. Ghelardini, C., Galeotti, N., Gualtieri, F., *et al.* The novel nootropic compound DM232 (unifiram) ameliorates memory impairment in mice and rats. *Drug Develop. Res.* **56**(1), 23-32 (2002).

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, 11/18/2022

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