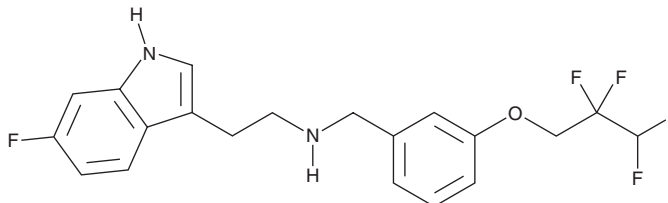


PRODUCT INFORMATION



Lu AE58054
Item No. 27335

CAS Registry No.: 467459-31-0
Formal Name: 6-fluoro-N-[[3-(2,2,3,3-tetrafluoropropoxy)phenyl]methyl]-1H-indole-3-ethanamine
MF: C₂₀H₁₉F₅N₂O
FW: 398.4
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 279 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

Lu AE58054 is supplied as a solid. A stock solution may be made by dissolving the Lu AE58054 in the solvent of choice, which should be purged with an inert gas. Lu AE58054 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of Lu AE58054 in these solvents is approximately 30 mg/ml.

Description

Lu AE58054 is an antagonist of the serotonin (5-HT) receptor subtype 5-HT₆ (K_i = 0.83 nM).¹ It is selective for 5-HT₆ over other 5-HT receptor subtypes (K_is = 250 to >10,000 nM) and 70 other targets but does bind to α_{1A}- and α_{1B}-adrenergic receptors (α_{1B}-ARs; K_is = 21 and 22 nM, respectively). Lu AE58054 (5-20 mg/kg) reverses phencyclidine-induced deficits in novel object recognition memory in rats. When administered in combination with the cholinesterase inhibitor rivastigmine (Item No. 14270), Lu AE58054 has an additive effect on the decreased number of slips and falls made by rats with bilateral striatal-dopaminergic and cortical-cholinergic system lesions, a model of Parkinson's disease motor disruption, in the Michigan complex motor control task.²

References

1. Arnt, J., Bang-Andersen, B., Grayson, B., *et al.* Lu AE58054, a 5-HT₆ antagonist, reverses cognitive impairment induced by subchronic phencyclidine in a novel object recognition test in rats. *Int. J. Neuropsychopharmacol.* **13**(8), 1021-1033 (2010).
2. Cherian, A.K., Kucinski, A., Wu, R., *et al.* Co-treatment with rivastigmine and idalopirdine reduces the propensity for falls in a rat model of falls in Parkinson's disease. *Psychopharmacology (Berl)* **236**(6), 1701-1715 (2019).

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, 12/15/2022

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