

PRODUCT INFORMATION



L-694,247 (trifluoroacetate salt)

Item No. 40639

Formal Name: N-[4-[[5-[3-(2-aminoethyl)-1H-indol-5-yl]-1,2,4-oxadiazol-3-yl]methyl]phenyl]-methanesulfonamide, trifluoroacetate salt

MF: C₂₀H₂₁N₅O₃S • XCF₃COOH

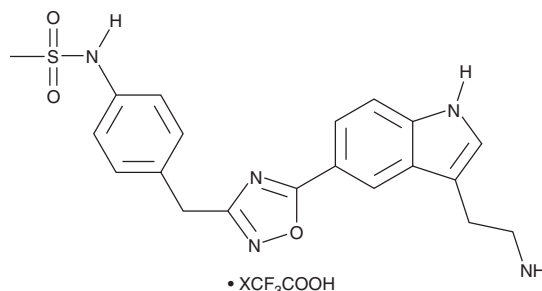
FW: 411.5

Purity: ≥95%

Supplied as: A crystalline 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

L-694,247 (trifluoroacetate salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the L-694,247 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. L-694,247 (trifluoroacetate salt) is slightly soluble (0.1-1 mg/ml) in acetonitrile.

L-694,247 (trifluoroacetate salt) is sparingly soluble (1-10 mg/ml) in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

L-694,247 is an agonist of the serotonin (5-HT) receptor subtype 5-HT_{1D}.¹ It is selective for 5-HT_{1D} over the 5-HT_{1A}, 5-HT_{1C}, 5-HT_{1E}, 5-HT₂, and 5-HT₃ receptors (K_is = 0.093, 2.29, 380, 2,187, 316, and >10,000 nM, respectively). L-694,247 inhibits adenylyl cyclase activity induced by forskolin (Item No. 11018) in guinea pig substantia nigra homogenates and potassium-induced 5-HT release in isolated guinea pig frontal cortex strips (EC₅₀s = 0.79 and 0.4 nM, respectively). It decreases water intake in dehydrated rats as well as prevents carbachol-, angiotensin II-, or isoproterenol-induced increases in water intake in normohydrated rats in a dose-dependent manner.² Intra-atrial administration of L-694,247 (0.00000125-0.1 µg/kg) decreases renal perfusion pressure but does not affect systolic blood pressure or heart rate in *in situ* autoperfused rat kidney infused with phenylephrine (Item Nos. 17205 | 18619).³

References

1. Beer, M.S., Stanton, J.A., Bevan, Y., *et al*. L-694,247: A potent 5-HT_{1D} receptor agonist. *Br. J. Pharmacol.* **110**(3), 1196-1200 (1993).
2. De Castro-e-Silva, E., Sarmento, C., Nascimento, T.A., *et al*. Effect of third ventricle administration of L-694,247, a selective 5-HT_{1D} receptor agonist, on water intake in rats. *Pharmacol. Biochem. Behav.* **57**(4), 749-754 (1997).
3. García-Pedraza, J.Á., García, M., Martín, M.L., *et al*. Pharmacological evidence that 5-HT_{1D} activation induces renal vasodilation by NO pathway in rats. *Clin. Exp. Pharmacol. Physiol.* **42**(6), 640-647 (2015).

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, 05/01/2024

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