

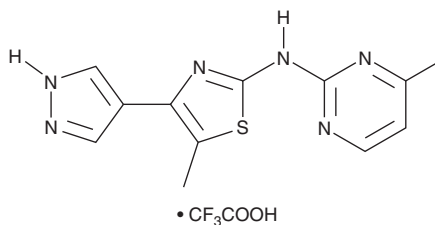
# PRODUCT INFORMATION



## ADX-88178 (trifluoroacetate salt)

Item No. 43178

**Formal Name:** 5-methyl-N-(4-methylpyrimidin-2-yl)-4-(1H-pyrazol-4-yl)thiazol-2-amine, trifluoroacetate salt  
**MF:** C<sub>12</sub>H<sub>12</sub>N<sub>6</sub>S • CF<sub>3</sub>COOH  
**FW:** 386.4  
**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

ADX-88178 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the ADX-88178 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. ADX-88178 (trifluoroacetate salt) is sparingly soluble (1-10 mg/ml) in DMSO.

### Description

ADX-88178 is a positive allosteric modulator (PAM) of metabotropic glutamate receptor 4 (mGluR4).<sup>1</sup> It enhances glutamate-induced intracellular calcium mobilization in HEK293 cells expressing human or rat mGluR4 (EC<sub>50</sub>s = 3.5 and 9.1 nM, respectively). ADX-88178 is selective for mGluR4 over mGluR1, mGluR2, mGluR3, mGluR5, mGluR7, and the GABA<sub>B</sub> receptor at 30 μM, mGluR6 at 10 μM, and mGluR8 (EC<sub>50</sub> = 2.2 μM). It is also selective for mGluR4 over a panel of 71 receptors, transporters, enzymes, and ion channels at 10 μM. ADX-88178 (10 mg/kg) reverses catalepsy induced by haloperidol (Item No. 12014) in rats. It enhances L-DOPA-induced decreases in motor impairments in 6-OHDA-lesioned rats in a dose-dependent manner. However, it does not reverse L-DOPA- and benserazide-induced dyskinesia in 6-OHDA-lesioned rats.<sup>2</sup>

### References

1. Le Poul, E., Boléa, C., Girard, F., *et al.* A potent and selective metabotropic glutamate receptor 4 positive allosteric modulator improves movement in rodent models of Parkinson's disease. *J. Pharmacol. Exp. Ther.* **343**(1), 167-177 (2012).
2. Finlay, C.J., Jackson, M.J., Fisher, R., *et al.* Metabotropic glutamate receptor 4 (mGlu4) positive allosteric modulators lack efficacy in rat and marmoset models of L-DOPA-induced dyskinesia. *J. Parkinsons Dis.* **14**(2), 245-259 (2024).

#### 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, 06/27/2025

#### 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