

PRODUCT INFORMATION



HTMT (maleate)

Item No. 29018

CAS Registry No.: 195867-54-0

Formal Name: 6-[[2-(1H-imidazol-5-yl)ethyl]amino]-N-[4-(trifluoromethyl)phenyl]-heptanamide, 2Z-butenedioate

Synonym: Histamine trifluoromethyl toluidine

MF: $C_{19}H_{25}F_3N_4O \cdot 2C_4H_4O_4$

FW: 614.6

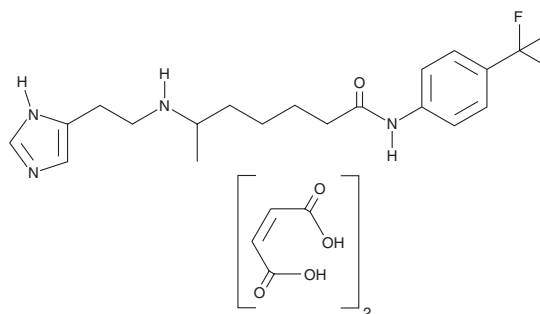
Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 250 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

HTMT (maleate) is supplied as a solid. A stock solution may be made by dissolving the HTMT (maleate) in the solvent of choice, which should be purged with an inert gas. HTMT (maleate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of HTMT (maleate) in these solvents is approximately 10, 25, and 30 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of HTMT (maleate) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of HTMT (maleate) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

HTMT is a histamine H_1 receptor agonist.¹ It also binds to the histamine H_4 receptor ($K_i = 1.2 \mu\text{M}$) but does not increase calcium mobilization in HEK293 cells expressing the receptor at $10 \mu\text{M}$.² HTMT increases intracellular calcium and inositol phosphate levels (EC_{50} s = 19 and $30 \mu\text{M}$, respectively) in human peripheral blood lymphocytes.³ It increases proliferation of immortalized mouse small, but not large, cholangiocytes, when used at a concentration of $10 \mu\text{M}$.⁴ HTMT ($10 \mu\text{g/kg}$ twice per day) increases levels of sheep red blood cell-induced IgG and IgM antibodies in rabbit serum.⁵ It induces scratching in mice when administered at an intradermal dose of $0.2 \mu\text{mol}$, an effect that is reduced by the H_1 receptor antagonist terfenadine (Item No. 20305).⁶

References

1. Shayo, C., Fernandez, N., Legnazzi, B.L., et al. *Mol. Pharmacol.* **60**(5), 1049-1056 (2001).
2. Zhu, Y., Michalovich, D., Wu, H., et al. *Mol. Pharmacol.* **59**(3), 434-441 (2001).
3. Qiu, R., Melmon, K.L., and Kahn, M.M. *J. Pharmacol. Exp. Ther.* **253**(3), 1245-1252 (1990).
4. Francis, H., Glaser, S., Demorrow, S., et al. *Am. J. Physiol. Cell Physiol.* **295**(2), C499-C513 (2008).
5. Tripathi, T., Shahid, M., Khan, H.M., et al. *Pharmacol. Rep.* **62**(5) 917-925 (2010).
6. Bell, J.K., McQueen, D.S., and Rees, J.L. *Br. J. Pharmacol.* **142**(2), 374-380 (2004).

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, 11/03/2022

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