

PRODUCT INFORMATION



Tegaserod (maleate)

Item No. 14692

CAS Registry No.: 189188-57-6

Formal Name: 2-[(5-methoxy-1H-indol-3-yl)methylene]-
N-pentyl-hydrazinecarboximidamide,
2Z-butenedioate

Synonyms: SDZ-HTF 919, Zelnorm

MF: $C_{16}H_{23}N_5O \cdot C_4H_4O_4$

FW: 417.5

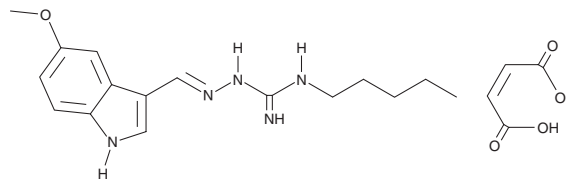
Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 279, 316 nm

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

Tegaserod (maleate) is supplied as a crystalline solid. A stock solution may be made by dissolving the tegaserod (maleate) in the solvent of choice. Tegaserod (maleate) is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of tegaserod (maleate) in these solvents is approximately 20 and 25 mg/ml, respectively.

Tegaserod (maleate) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, tegaserod (maleate) should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Tegaserod (maleate) has a solubility of approximately 0.03 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Tegaserod (maleate) is a potent agonist of the serotonin 4 receptor (5-HT_4 , $K_i = 3.9\text{-}31$ nM).^{1,2} As 5-HT_4 receptors are associated with cholinergic nerves lining the colon, tegaserod (maleate) stimulates colonic motility and transit *in vivo*, in both healthy and diseased gastrointestinal tracts.^{3,4} However, it lacks selectivity, since it also acts as an antagonist for the 5-HT_{2A} , 5-HT_{2B} , and 5-HT_{2C} receptors ($K_i = 32, 3.9,$ and 100 nM, respectively).⁴⁻⁶ It is a weak agonist of 5-HT_3 ($K_i \leq 10$ μM).⁷

References

1. Long, D.D., Armstrong, S.R., Beattie, D.T., *et al.* *Bioorg. Med. Chem. Lett.* **22**(19), 6048-6052 (2012).
2. Hedner, E., Sjögren, M., Frändberg, P.-A., *et al.* *J. Nat. Prod.* **69**(10), 1421-1424 (2006).
3. Nguyen, A., Camilleri, M., Kost, L.J., *et al.* *J. Pharmacol. Exp. Ther.* **280**(3), 1270-1276 (1997).
4. Tack, J., Camilleri, M., Chang, L., *et al.* *Aliment. Pharmacol. Ther.* **35**(7), 745-767 (2012).
5. Beattie, D.T., Smith, J.A.M., Marquess, D., *et al.* *Br. J. Pharmacol.* **143**(5), 549-560 (2004).
6. McCullough, J.L., Armstrong, S.R., Hegde, S.S., *et al.* *Pharmacol. Res.* **53**(4), 353-358 (2006).
7. Lemaître, S., Lepailleur, A., Bureau, R., *et al.* *Bioorg. Med. Chem.* **17**(6), 2607-2622 (2009).

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, 11/23/2022

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