

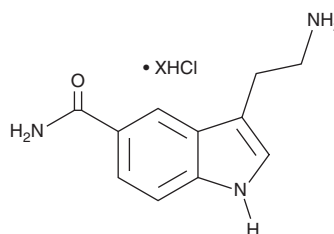
# PRODUCT INFORMATION



## 5-Carboxamidotryptamine (hydrochloride)

Item No. 45295

**CAS Registry No.:** 467451-77-0  
**Formal Name:** 3-(2-aminoethyl)-1H-indole-5-carboxamide, hydrochloride  
**Synonyms:** 5-CT  
**MF:** C<sub>11</sub>H<sub>13</sub>N<sub>3</sub>O • XHCl  
**FW:** 203.2  
**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

5-Carboxamidotryptamine (hydrochloride) is supplied as a solid. Aqueous solutions of 5-carboxamidotryptamine (hydrochloride) can be prepared by directly dissolving the solid in aqueous buffers. 5-Carboxamidotryptamine (hydrochloride) is soluble (≥10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

### Description

5-Carboxamidotryptamine is an agonist of the serotonin (5-HT) receptor subtypes 5-HT<sub>1</sub> and 5-HT<sub>7</sub>.<sup>1-3</sup> It binds to the 5-HT<sub>1D</sub> receptor in postmortem human cortical tissue (IC<sub>50</sub>s = 1.82 and 2,041 nM for the high- and low-affinity binding sites, respectively) and the 5-HT<sub>7</sub> receptor in HEK293 cells expressing 5-HT<sub>7</sub> (K<sub>i</sub> = 0.79 nM) in radioligand binding assays.<sup>3,4</sup> 5-Carboxamidotryptamine induces hypotension and bradycardia and increases hindlimb vascular conductance in anesthetized rats when administered via injection into the rostral ventrolateral medulla.<sup>5</sup>

### References

1. Hoyer, D. and Schoeffter, P. 5-HT receptors: Subtypes and second messengers. *J. Recept. Res.* **11(1-4)**, 197-214 (1991).
2. Zifa, E. and Fillion, G. 5-Hydroxytryptamine receptors. *Pharm. Rev.* **44(3)**, 401-458 (1992).
3. Wood, M., Chaubey, M., Atkinson, P., et al. Antagonist activity of meta-chlorophenylpiperazine and partial agonist activity of 8-OH-DPAT at the 5-HT<sub>7</sub> receptor. *Eur. J. Pharm.* **396(1)**, 1-8 (2000).
4. Beer, M.S., Stanton, J.A., Bevan, Y., et al. An investigation of the 5-HT<sub>1D</sub> receptor binding affinity of 5-hydroxytryptamine, 5-carboxamidotryptamine and sumatriptan in the central nervous system of seven species. *Eur. J. Pharm.* **213(2)**, 193-197 (1992).
5. Lovick, T.A. Systemic and regional haemodynamic responses to microinjection of 5-HT agonists in the rostral ventrolateral medulla in the rat. *Neurosci. Lett.* **107(1-3)**, 157-161 (1989).

#### 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, 05/06/2026

#### 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