

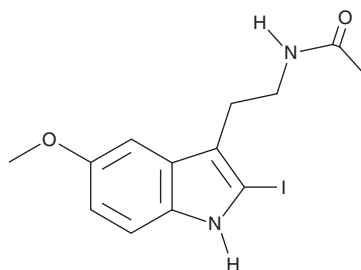
# PRODUCT INFORMATION



## 2-Iodomelatonin

Item No. 19711

**CAS Registry No.:** 93515-00-5  
**Formal Name:** N-[2-(2-iodo-5-methoxy-1H-indol-3-yl)ethyl]-acetamide  
**MF:** C<sub>13</sub>H<sub>15</sub>IN<sub>2</sub>O<sub>2</sub>  
**FW:** 358.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 214, 279 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

2-Iodomelatonin is supplied as a crystalline solid. A stock solution may be made by dissolving the 2-iodomelatonin in the solvent of choice, which should be purged with an inert gas. 2-Iodomelatonin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 2-iodomelatonin in ethanol is approximately 20 mg/ml and 30 mg/ml in DMSO and DMF.

2-Iodomelatonin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 2-iodomelatonin should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 2-Iodomelatonin has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

2-Iodomelatonin is a potent agonist of melatonin receptor 1 (MT<sub>1</sub>; K<sub>i</sub> = 28 pM) that is 5-fold selective for MT<sub>1</sub> over MT<sub>2</sub>.<sup>1,2</sup> It inhibits forskolin-stimulated cAMP production in CHO cells expressing human MT<sub>1</sub> 30-fold more potently than melatonin (Item No. 14427) with an EC<sub>50</sub> value of 11 pM. 2-Iodomelatonin has been used to characterize the role of MT<sub>1</sub> in melatonin-mediated signaling.<sup>3,4</sup>

### References

1. Dubocovich, M.L., Masana, M.I., Iacob, S., *et al.* Melatonin receptor antagonists that differentiate between the human Mel<sub>1a</sub> and Mel<sub>1b</sub> recombinant subtypes are used to assess the pharmacological profile of the rabbit retina ML<sub>1</sub> presynaptic heteroreceptor. *Naunyn-Schmiedeberg's Arch. Pharmacol.* **355**(3), 365-375 (1997).
2. Browning, C., Beresford, I., Fraser, N., *et al.* Pharmacological characterization of human recombinant melatonin mt<sub>1</sub> and MT<sub>2</sub> receptors. *Br. J. Pharmacol.* **129**(5), 877-886 (2000).
3. Cecon, E., Chen, M., Marçola, M., *et al.* Amyloid β peptide directly impairs pineal gland melatonin synthesis and melatonin receptor signaling through the ERK pathway. *FASEB J.* **29**(6), 2566-2582 (2015).
4. Adamah-Biassi, E.B., Zhang, Y., Jung, H., *et al.* Distribution of MT<sub>1</sub> melatonin receptor promoter-driven RFP expression in the brains of BAC C3H/HeN transgenic mice. *J. Histochem. Cytochem.* **62**(1), 70-84 (2014).

#### 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, 02/06/2025

#### 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