

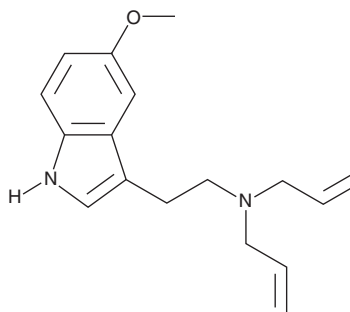
# PRODUCT INFORMATION



## 5-methoxy DALT

Item No. 10729

**CAS Registry No.:** 928822-98-4  
**Formal Name:** 5-methoxy-N,N-di-2-propen-1-yl-1H-indole-3-ethanamine  
**Synonyms:** N,N-Diallyl-5-Methoxytryptamine  
**MF:** C<sub>17</sub>H<sub>22</sub>N<sub>2</sub>O  
**FW:** 270.4  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 224, 279 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

5-methoxy DALT is supplied as a crystalline solid. A stock solution may be made by dissolving the 5-methoxy DALT in the solvent of choice, which should be purged with an inert gas. 5-methoxy DALT is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide (DMF). The solubility of 5-methoxy DALT in these solvents is approximately 1.5, 1, 15, and 20 mg/ml, respectively.

5-methoxy DALT is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 5-methoxy DALT should first be dissolved in DMF and then diluted with the aqueous buffer of choice. 5-methoxy DALT has a solubility of approximately 0.15 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

5-methoxy DALT is a tryptamine derivative recently used as a component in 'bath salts' that act directly on monoamine receptors (similar to 3,4-methylenedioxyamphetamine) to produce its psychoactive effects.<sup>1</sup> The physiological, neurological, and toxicological actions of 5-methoxy DALT have not been fully characterized. This compound is intended as an analytical standard for the forensic analysis of samples that may contain this compound.

### Reference

1. Nonaka, R., Nagai, F., Ogata, A., *et al.* *In vitro* screening of psychoactive drugs by [<sup>35</sup>S]GTPγS binding in rat brain membranes. *Biol. Pharm. Bull.* **30(12)**, 2328-33 (2007).

#### 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, 05/17/2022

#### 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