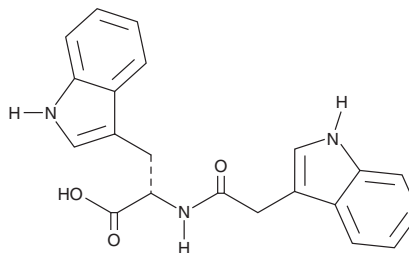


# PRODUCT INFORMATION

## (2-(1H-Indol-3-yl)acetyl)-L-Tryptophan

Item No. 43104

**CAS Registry No.:** 57105-53-0  
**Synonyms:** (2-(1H-Indol-3-yl)acetyl)-L-Trp, IAA-Trp, IATrp, Indole-3-acetyl-L-tryptophan  
**MF:** C<sub>21</sub>H<sub>19</sub>N<sub>3</sub>O<sub>3</sub>  
**FW:** 361.4  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Synthetic



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

### Laboratory Procedures

(2-(1H-Indol-3-yl)acetyl)-L-Tryptophan is supplied as a solid. A stock solution may be made by dissolving the (2-(1H-Indol-3-yl)acetyl)-L-tryptophan in the solvent of choice, which should be purged with an inert gas. (2-(1H-Indol-3-yl)acetyl)-L-Tryptophan is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

### Description

(2-(1H-Indol-3-yl)acetyl)-L-tryptophan is an active metabolite of the auxin plant growth regulator indole-3-acetic acid (IAA; Item No. 16954).<sup>1,2</sup> It induces coleoptile elongation in *A. sativa* seedlings ( $EC_{50} = 6.5 \mu M$ ).<sup>2</sup> (2-(1H-Indol-3-yl)acetyl)-L-tryptophan inhibits IAA-induced inhibition of *A. thaliana* root growth in a concentration-dependent manner.<sup>1</sup>

### References

1. Staswick, P.E. The tryptophan conjugates of jasmonic and indole-3-acetic acids are endogenous auxin inhibitors. *Plant. Physiol.* **150**(3), 1310-1321 (2009).
2. Feung, C.-S., Hamilton, R.H., and Mumma, R.O. Metabolism of indole-3-acetic acid. *Plant. Physiol.* **59**(1), 91-93 (1977).

#### 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, 04/11/2025

#### 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