

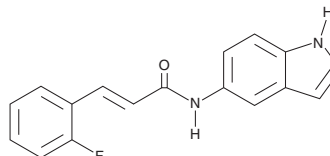
# PRODUCT INFORMATION



## (E)-3-(2-Fluorophenyl)-N-(1H-indol-5-yl)acrylamide

Item No. 43022

CAS Registry No.: 3094045-97-0  
Formal Name: 3-(2-fluorophenyl)-N-1H-indol-5-yl-  
(2E)-propenamide  
MF: C<sub>17</sub>H<sub>13</sub>FN<sub>2</sub>O  
FW: 280.3  
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

(E)-3-(2-Fluorophenyl)-N-(1H-indol-5-yl)acrylamide is supplied as a solid. A stock solution may be made by dissolving the (E)-3-(2-fluorophenyl)-N-(1H-indol-5-yl)acrylamide in the solvent of choice, which should be purged with an inert gas. (E)-3-(2-Fluorophenyl)-N-(1H-indol-5-yl)acrylamide is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

### Description

(E)-3-(2-Fluorophenyl)-N-(1H-indol-5-yl)acrylamide is an inhibitor of monoamine oxidase B (MAO-B) and an inducer of nuclear factor erythroid 2-related factor 2 (Nrf2).<sup>1</sup> It inhibits MAO-B (IC<sub>50</sub> = 3.61 μM) and increases levels of Nrf2 in a reporter assay using AREc32 cells (IC<sub>50</sub> = 0.361 μM). (E)-3-(2-Fluorophenyl)-N-(1H-indol-5-yl)acrylamide scavenges peroxy radicals in an oxygen radical absorbance capacity (ORAC) assay (IC<sub>50</sub> = 2.77 μM) and inhibits LPS-induced increases in nitrite levels in BV-2 microglial cells (EC<sub>50</sub> = 11.6 μM). It inhibits 6-OHDA- or rotenone- and oligomycin A-induced decreases in viability in SH-SY5Y neuroblastoma cells when used at a concentration of 1 μM.

### Reference

1. Duarte, P., Sanchez-Porro, F.J., Crisman, E., *et al.* Network-based drug optimization toward the treatment of Parkinson's disease: NRF2, MAO-B, oxidative stress, and chronic neuroinflammation. *J. Med. Chem.* **68**(3), 3495-3517 (2025).

#### 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, 02/03/2026

#### 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