

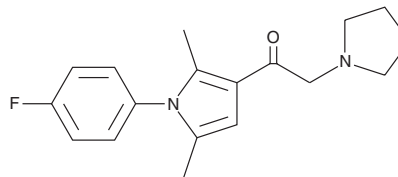
# PRODUCT INFORMATION



## IU1

Item No. 10617

**CAS Registry No.:** 314245-33-5  
**Formal Name:** 1-[1-(4-fluorophenyl)-2,5-dimethyl-1H-pyrrol-3-yl]-2-(1-pyrrolidinyl)-ethanone  
**MF:** C<sub>18</sub>H<sub>21</sub>FN<sub>2</sub>O  
**FW:** 300.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 247, 289 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

IU1 is supplied as a crystalline solid. A stock solution may be made by dissolving the IU1 in the solvent of choice. IU1 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of IU1 in these solvents is approximately 20 mg/ml.

IU1 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, IU1 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. IU1 has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Polyubiquitinylation designates proteins for proteasomal degradation. USP14 is one of several deubiquitinating enzymes (DUBs) responsible for shortening proteasome-bound ubiquitin chains, which is known to antagonize the degradation of ubiquitin-protein conjugates.<sup>1</sup> IU1 is a reversible, small molecule inhibitor of deubiquitination by USP14 that demonstrates an IC<sub>50</sub> value of 4-5 μM. IU1 selectively stimulates ubiquitin-dependent protein degradation *in vitro* at 34 μM and in MEF cells at 50 μM. At 75 μM, IU1 reduces accumulation of oxidized proteins in HEK293 cells, alleviating cytotoxicity induced by oxidative stress.<sup>1</sup>

### Reference

1. Lee, B.-H., Lee, M.J., Park, S., *et al.* Enhancement of proteasome activity by a small-molecule inhibitor of USP14. *Nature* **467**, 179-184 (2010).

#### 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, 11/08/2022

#### 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