

IU1 (1-[1-(4-Fluorophenyl)-2,5-dimethyl-1H-pyrrol-3-yl]-2-(1-pyrrolidinyl)ethanone

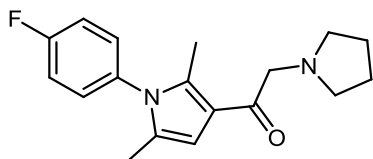
Cat. # SI9629

Background: IU1 is a cell-permeable, reversible and selective inhibitor of proteasome-associated Usp14 ($IC_{50} < 4 \mu\text{M}$). Enhances proteasome activity and substrate proteolysis. IU1 promotes cell survival under stress conditions and selectively stimulates ubiquitin-dependent protein degradation *in vitro*.

Application: Inhibition of USP14.

Product Information

| | |
|--------------------------|--|
| Purity: | > 98% (TLC); NMR (Conforms) |
| Molecular Weight: | 300.4 Da |
| Physical State: | Beige powder |
| Quantity: | 10 mg |
| Solubility: | DMSO (30mg/ml) |
| Storage: | Store desiccated as supplied at 4°C for up to 2 years. Store solutions at -20°C for up to one month. Protect from light. |



Formula: C₁₈H₂₁FN₂O

References

1. Lee et al (2010) Enhancement of proteasome activity by a small-molecule inhibitor of USP14. Nature 467 179. PMID: 20829789.
2. Lee et al (2010) Trimming of ubiquitin chains by proteasome-associated deubiquitinating enzymes. Mol.Cell.Proteomics 10 R110.

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses
Copyright © 2011 LifeSensors, Inc. All Rights Reserved