

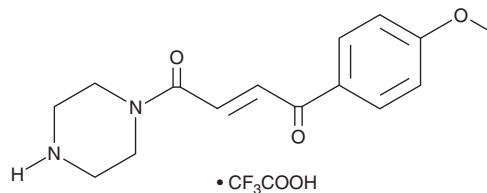
# PRODUCT INFORMATION



**JP-2-196**

Item No. 42342

**CAS Registry No.:** 2913587-31-0  
**Formal Name:** 1-(4-methoxyphenyl)-4-(1-piperazinyl)-2-butene-1,4-dione  
2,2,2-trifluoroacetate  
**MF:** C<sub>15</sub>H<sub>18</sub>N<sub>2</sub>O<sub>3</sub> • CF<sub>3</sub>COOH  
**FW:** 388.3  
**Purity:** ≥90%  
**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

JP-2-196 is supplied as a solid. A stock solution may be made by dissolving the JP-2-196 in the solvent of choice, which should be purged with an inert gas. JP-2-196 is sparingly (1-10 mg/ml) soluble in DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of JP-2-196 can be prepared by directly dissolving the solid in aqueous buffers. JP-2-196 is sparingly (1-10 mg/ml) soluble in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

## Description

JP-2-196 is a molecular glue handle.<sup>1</sup> It binds to the E3 ubiquitin ligases RNF126, LRSAM1, RNF40, MID2, RNF219, and RNF14 in an immunoprecipitation (IP) assay when used at a concentration of 10 μM. JP-2-196 has been used in the synthesis of various molecular glue degraders.

## Reference

1. Toriki, E.S., Papatzimas, J.W., Nishikawa, K., *et al.* Rational chemical design of molecular glue degraders. *ACS Cent. Sci.* **9**(5), 915-926 (2023).

### 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, 01/27/2025

## 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