

# PRODUCT INFORMATION



## STING PROTAC SP23

Item No. 39274

**CAS Registry No.:** 2762552-74-7  
**Formal Name:** (2E)-N1-[6-[[2-(2,6-dioxo-3-piperidinyl)-2,3-dihydro-1,3-dioxo-1H-isoindol-4-yl]amino]hexyl]-N4-[4-[[[(5-nitro-2-furanyl)carbonyl]amino]phenyl]-2-butenediamide

**Synonyms:** PROTAC Stimulator of Interferon Genes Degradar-1, PROTAC STING Degradar-1, Proteolysis-targeting Chimera Stimulator of Interferon Genes Degradar-1, Proteolysis-targeting Chimera STING Degradar-1, Stimulator of Interferon Genes PROTAC SP23, Stimulator of Interferon Genes Proteolysis-targeting Chimera SP23, STING Proteolysis-targeting Chimera SP23

**MF:** C<sub>34</sub>H<sub>33</sub>N<sub>7</sub>O<sub>10</sub>

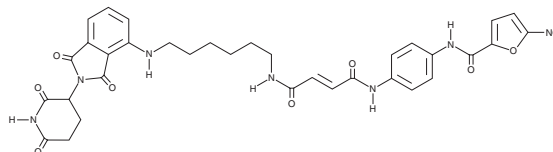
**FW:** 699.7

**Purity:** ≥95%

**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

STING PROTAC SP23 is supplied as a solid. A stock solution may be made by dissolving the STING PROTAC SP23 in the solvent of choice, which should be purged with an inert gas. STING PROTAC SP23 is soluble in acetonitrile.

### Description

STING PROTAC SP23 is a proteolysis-targeting chimera (PROTAC) containing the stimulator of interferon genes (STING) inhibitor C-170 (Item No. 30157) conjugated to the cereblon (CRBN) inhibitor pomalidomide (Item No. 19877) via an  $\alpha,\beta$ -unsaturated carbonyl linker.<sup>1</sup> It degrades STING in THP-1 cells with a half-maximal degradation concentration (DC<sub>50</sub>) of 3.2  $\mu$ M, an effect that can be blocked by the proteasome inhibitor MG132 but not the lysosome inhibitor bafilomycin. STING PROTAC SP23 (2.5-20  $\mu$ M) decreases cGAMP-induced increases in the levels of IFN- $\beta$ , IL-6, and chemokine (C-X-C motif) ligand 10 (CXCL10) in THP-1 cells. It also reduces kidney weight, improves kidney function, and increases the survival rate in a mouse model of acute kidney injury induced by the DNA-crosslinking agent cisplatin (Item No. 13119) when administered at doses of 30 and 60 mg/kg.

### Reference

1. Liu, J., Yuan, L., Ruan, Y., *et al.* Novel CRBN-recruiting proteolysis-targeting chimeras as degraders of stimulator of interferon genes with in vivo anti-inflammatory efficacy. *J. Med. Chem.* **65(9)**, 6593-6611 (2022).

#### 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, 02/13/2024

#### 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