

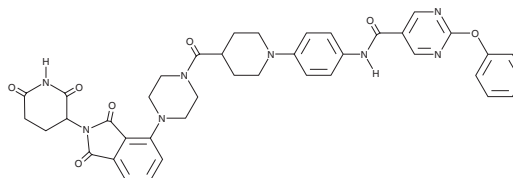
# PRODUCT INFORMATION



## PROTAC(H-PGDS)-7

Item No. 35561

**CAS Registry No.:** 2761281-50-7  
**Formal Name:** N-[4-[4-[[4-[2-(2,6-dioxo-3-piperidiny)]-2,3-dihydro-1,3-dioxo-1H-isoindol-4-yl]-1-piperazinyl] carbonyl]-1-piperidiny] phenyl]-2-phenoxy-5-pyrimidinecarboxamide  
**Synonym:** Proteolysis-targeting Chimera (Prostaglandin D Synthase (hematopoietic-type))-7



**MF:** C<sub>40</sub>H<sub>38</sub>N<sub>8</sub>O<sub>7</sub>  
**FW:** 742.8  
**Purity:** ≥95%  
**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

PROTAC(H-PGDS)-7 is supplied as a crystalline solid. A stock solution may be made by dissolving the PROTAC(H-PGDS)-7 in the solvent of choice, which should be purged with an inert gas. PROTAC(H-PGDS)-7 is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

### Description

PROTAC(H-PGDS)-7 is a proteolysis-targeting chimera (PROTAC) comprised of the hematopoietic prostaglandin D synthase (H-PGDS) inhibitor TFC-007 (Item No. 17971) and the cereblon inhibitor pomalidomide (Item No. 19877).<sup>1</sup> It induces degradation of H-PGDS in KU812 cells with a 50% degradation concentration (DC<sub>50</sub>) value of 17.3 pM, as well as decreases the production of prostaglandin D<sub>2</sub> (PGD<sub>2</sub>) in the same cells.

### Reference

1. Yokoo, H., Shibata, N., Endo, A., *et al.* Discovery of a highly potent and selective degrader targeting hematopoietic prostaglandin D synthase via in silico design. *J. Med. Chem.* **64**(21), 15868-15882 (2021).

#### 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, 10/09/2025

#### 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