

# PRODUCT INFORMATION

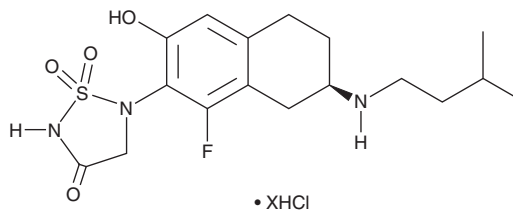


## ABBV-CLS-484 (hydrochloride)

Item No. 40169

**Formal Name:** 5-[(7R)-1-fluoro-5,6,7,8-tetrahydro-3-hydroxy-7-[(3-methylbutyl)amino]-2-naphthalenyl]-1,2,5-thiadiazolidin-3-one, 1,1-dioxide, hydrochloride

**Synonym:** Osunprotafib  
**MF:** C<sub>17</sub>H<sub>24</sub>FN<sub>3</sub>O<sub>4</sub>S • XHCl  
**FW:** 385.5  
**Purity:** ≥98%  
**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

ABBV-CLS-484 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the ABBV-CLS-484 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. ABBV-CLS-484 (hydrochloride) is slightly soluble in acetonitrile.

ABBV-CLS-484 (hydrochloride) is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

### Description

ABBV-CLS-484 is a dual inhibitor of protein tyrosine phosphatase 1N (PTP1N) and PTP2N (IC<sub>50</sub>s = 2.5 and 1.8 nM, respectively).<sup>1</sup> It is selective for PTP1N and PTP2N over PTP6N and PTP11N (IC<sub>50</sub>s = >10 μM for both) but does inhibit PTP9N (IC<sub>50</sub> = 15 nM). ABBV-CLS-484 (1 μM) potentiates IFN-γ-induced decreases in the proliferation of B16-F1 mouse melanoma cells. *In vivo*, ABBV-CLS-484 (20 mg/kg per day), in combination with an anti-PD-L1 antibody, reduces tumor volume in a B16-F1 melanoma murine model. ABBV-CLS-484 alone increases the number of CD8<sup>+</sup> effector T cells, natural killer (NK) T cells, neutrophils, and monocytes expressing high levels of IFN-γ-induced genes, as well as increases M1 and decreases M2 macrophages, in the tumor microenvironment in the same mouse model when administered at a dose of 20 mg/kg per day.

### Reference

1. Baumgartner, C.K., Ebrahimi-Nik, H., Iracheta-Vellve, A., *et al.* The PTPN2/PTPN1 inhibitor ABBV-CLS-484 unleashes potent anti-tumour immunity. *Nature* **622(7984)**, 850-862 (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, 01/29/2024

#### 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