

# PRODUCT INFORMATION

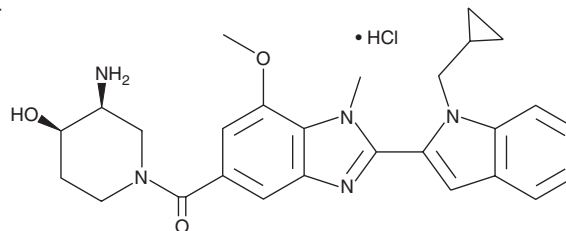


## GSK484 (hydrochloride)

Item No. 17488

**CAS Registry No.:** 1652591-81-5  
**Formal Name:** [(3S,4R)-3-amino-4-hydroxy-1-piperidinyl][2-[1-(cyclopropylmethyl)-1H-indol-2-yl]-7-methoxy-1-methyl-1H-benzimidazol-5-yl]-methanone, monohydrochloride

**MF:** C<sub>27</sub>H<sub>31</sub>N<sub>5</sub>O<sub>3</sub> • HCl  
**FW:** 510.0  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 227, 242, 304 nm  
**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

GSK484 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the GSK484 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. GSK484 (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). It is also soluble in water. The solubility of GSK484 (hydrochloride) in ethanol and DMSO is approximately 25 mg/ml, approximately 30 mg/ml in DMF, and approximately 10 mg/ml in water. We do not recommend storing the aqueous solution for more than one day.

### Description

GSK484 is a reversible inhibitor of PAD4 (IC<sub>50</sub> = 50 nM) that binds to the low-calcium form of the enzyme.<sup>1</sup> It is selective for PAD4 over PAD1-3. GSK484 blocks the citrullination of PAD4 target proteins in human neutrophils and inhibits the formation of neutrophil extracellular traps in both mouse and human neutrophils.<sup>1</sup> It exhibits favorable pharmacokinetic profiles in mouse and rat. See the Structural Genomics Consortium (SGC) website for more information.

### Reference

1. Lewis, H.D., Liddle, J., Coote, J.E., *et al.* Inhibition of PAD4 activity is sufficient to disrupt mouse and human NET formation. *Nat. Chem. Biol.* **11**(3), 189-191 (2015).

#### 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, 10/16/2023

#### 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](http://WWW.CAYMANCHEM.COM)