

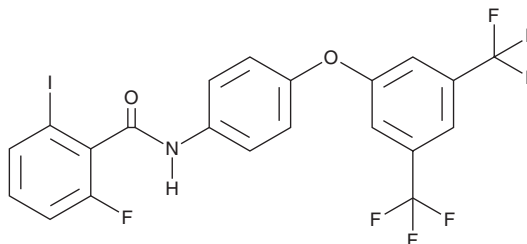
# PRODUCT INFORMATION



## CU-115

Item No. 29815

**CAS Registry No.:** 2471982-20-2  
**Formal Name:** N-[4-[3,5-bis(trifluoromethyl)phenoxy]phenyl]-2-fluoro-6-iodobenzamide  
**MF:** C<sub>21</sub>H<sub>11</sub>F<sub>7</sub>INO<sub>2</sub>  
**FW:** 569.2  
**Purity:** ≥98%  
**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

CU-115 is supplied as a crystalline solid. A stock solution may be made by dissolving the CU-115 in the solvent of choice, which should be purged with an inert gas. CU-115 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CU-115 in ethanol is approximately 0.25 mg/ml and approximately 10 mg/ml in DMSO and DMF.

CU-115 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CU-115 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CU-115 has a solubility of approximately 0.15 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

CU-115 is a toll-like receptor 8 (TLR8) antagonist (IC<sub>50</sub> = 1.04 μM).<sup>1</sup> It is selective for TLR8 over TLR7 (IC<sub>50</sub> = >50 μM). CU-115 inhibits increases in type I IFN transcriptional activity induced by the ssRNA nucleic acid ligands 3p-hpRNA or G3-YSD in a luciferase reporter assay when used at concentrations of 5 and 20 μM. It decreases TNF-α and IL-1β production induced by R-848 (Item No. 14806) in THP-1 cells when used at a concentration of 5 μM.

### Reference

1. Padilla-Salinas, R., Anderson, R., Sakaniwa, K., *et al.* Discovery of novel small molecule dual inhibitors targeting toll-like receptors 7 and 8. *J. Med. Chem.* **62**(22), 10221-10244 (2019).

#### 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, 03/20/2024

#### 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