

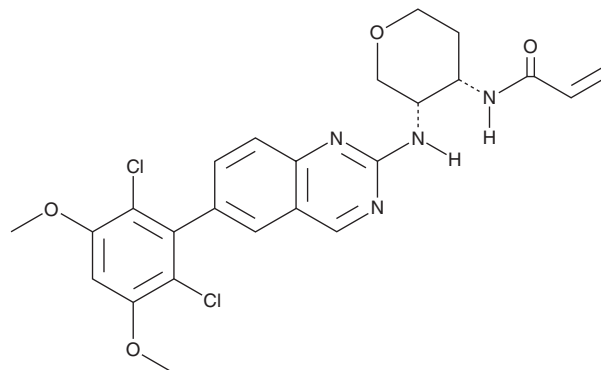
# PRODUCT INFORMATION



## BLU-554

Item No. 22438

**CAS Registry No.:** 1707289-21-1  
**Formal Name:** N-[(3S,4S)-3-[[6-(2,6-dichloro-3,5-dimethoxyphenyl)-2-quinazolinyl]amino]tetrahydro-2H-pyran-4-yl]-2-propenamide  
**MF:** C<sub>24</sub>H<sub>24</sub>Cl<sub>2</sub>N<sub>4</sub>O<sub>4</sub>  
**FW:** 503.4  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 246, 369 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

BLU-554 is supplied as a crystalline solid. A stock solution may be made by dissolving the BLU-554 in the solvent of choice. BLU-554 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of BLU-554 in ethanol is approximately 0.2 mg/ml and approximately 20 mg/ml in DMSO and DMF.

BLU-554 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BLU-554 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. BLU-554 has a solubility of approximately 0.3 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

BLU-554 is a potent inhibitor of recombinant fibroblast growth factor receptor 4 (FGFR4; IC<sub>50</sub> < 10 nM).<sup>1</sup> It also inhibits phosphorylation of ERK (IC<sub>50</sub> < 10 nM), a marker of FGFR4 inhibition, in MDA-MB-453 breast cancer cells.

### Reference

1. Bifulco, N., Jr., DiPietro, L.V., Hodous, B.L., *et al.* Inhibitors of the fibroblast growth factor receptor. United States, Blueprint Medicines Corp. (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, 10/26/2022

#### 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