

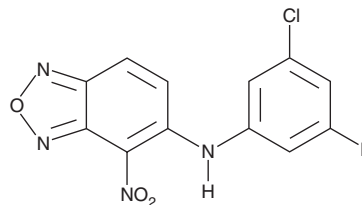
# PRODUCT INFORMATION



## HIF-2 $\alpha$ Antagonist 2

Item No. 16971

**CAS Registry No.:** 1422955-31-4  
**Formal Name:** N-(3-chloro-5-fluorophenyl)-4-nitro-2,1,3-benzoxadiazol-5-amine  
**MF:** C<sub>12</sub>H<sub>6</sub>ClFN<sub>4</sub>O<sub>3</sub>  
**FW:** 308.7  
**Purity:**  $\geq$ 98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

HIF-2 $\alpha$  antagonist 2 is supplied as a solid. A stock solution may be made by dissolving the HIF-2 $\alpha$  antagonist 2 in the solvent of choice, which should be purged with an inert gas. HIF-2 $\alpha$  antagonist 2 is soluble in the organic solvent DMSO at a concentration of approximately 30 mg/ml.

### Description

HIF-2 $\alpha$  antagonist 2 is an inhibitor of hypoxia-inducible factor-2 $\alpha$  (HIF-2 $\alpha$ ) PAS-B domain heterodimerization.<sup>1</sup> It binds to the HIF-2 $\alpha$  PAS-B domain ( $K_d = 81$  nM) to inhibit heterodimerization with the aryl hydrocarbon receptor nuclear translocator (ARNT) PAS-B domain in cell-free assays when used at concentrations ranging from 0.1 to 100  $\mu$ M. HIF-2 $\alpha$  antagonist 2 (0.1-30  $\mu$ M) decreases expression of the HIF-2 $\alpha$  target gene VEGFA in 786-O renal carcinoma cells.

### Reference

1. Scheuermann, T.H., Li, Q., Ma, H.-W., *et al.* Allosteric inhibition of hypoxia inducible factor-2 with small molecules. *Nat. Chem. Biol.* **9**(4), 271-276 (2013).

#### 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, 10/18/2022

#### 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