

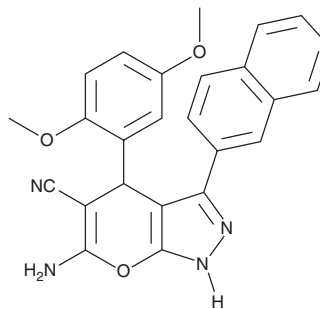
# PRODUCT INFORMATION



## RBC8

Item No. 21281

**CAS Registry No.:** 361185-42-4  
**Formal Name:** 6-amino-4-(2,5-dimethoxyphenyl)-1,4-dihydro-3-(2-naphthalenyl)-pyrano[2,3-c]pyrazole-5-carbonitrile  
**MF:** C<sub>25</sub>H<sub>20</sub>N<sub>4</sub>O<sub>3</sub>  
**FW:** 424.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 252 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

RBC8 is supplied as a crystalline solid. A stock solution may be made by dissolving the RBC8 in the solvent of choice, which should be purged with an inert gas. RBC8 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of RBC8 in ethanol is approximately 0.25 mg/ml and approximately 30 mg/ml in DMSO and DMF.

RBC8 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, RBC8 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. RBC8 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

RBC8 is a small molecule inhibitor of the Ras-like GTPases RalA and RalB that prevents binding of the GTPase to the Ral-binding protein.<sup>1</sup> RBC8 has been shown to prevent Ral-mediated spreading of murine embryonic fibroblasts, block anchorage-independent growth of H2122 and H358 lung cancer cell lines (IC<sub>50</sub>s = 3.5 and 3.4 μM, respectively), and inhibit the growth of H2122 lung cancer xenografts in mice (50 mg/kg i.p.).<sup>1</sup>

### Reference

1. Yan, C., Liu, D., Li, L., *et al.* Discovery and characterization of small molecules that target the GTPase Ral. *Nature* (2014).

#### 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, 12/01/2022

#### 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