

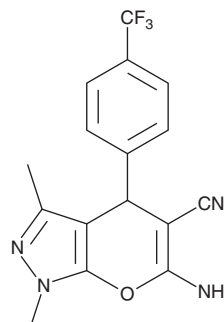
# PRODUCT INFORMATION



## BQU57

Item No. 16960

**CAS Registry No.:** 1637739-82-2  
**Formal Name:** 6-amino-1,4-dihydro-1,3-dimethyl-4-[4-(trifluoromethyl)phenyl]-pyrano[2,3-c]pyrazole-5-carbonitrile  
**MF:** C<sub>16</sub>H<sub>13</sub>F<sub>3</sub>N<sub>4</sub>O  
**FW:** 334.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 254 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

BQU57 is supplied as a crystalline solid. A stock solution may be made by dissolving the BQU57 in the solvent of choice, which should be purged with an inert gas. BQU57 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of BQU57 in ethanol is approximately 5 mg/ml and approximately 30 mg/ml in DMSO and DMF.

BQU57 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BQU57 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. BQU57 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

BQU57 is an inhibitor of the Ras-like GTPases RalA and RalB.<sup>1</sup> It binds to RalB when RalB is bound to GDP with a K<sub>d</sub> value of 7.7 μM, as determined by isothermal titration calorimetry (ITC). BQU57 (10 μM) decreases RalA and RalB activity in H2122 and H358 cells. It also inhibits colony formation of Ral-dependent H2122 and H358 cells (IC<sub>50</sub>s = 2 and 1.3 μM, respectively) but not Ral-independent H460 and Calu-6 cells. BQU57 (10, 20, and 50 mg/kg per day) decreases RalA and RalB, but not Ras or RhoA, activity and reduces tumor growth in a dose-dependent manner in an H2122 mouse xenograft model.

### Reference

1. Yan, C., Liu, D., Li, L., *et al.* Discovery and characterization of small molecules that target the GTPase Ral. *Nature* **515**(7527), 443-447 (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, 10/12/2022

#### 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