

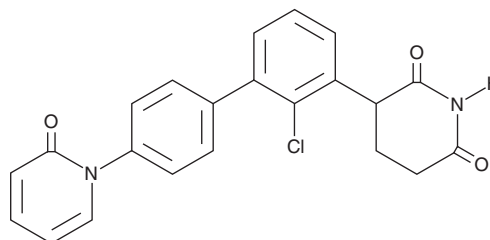
# PRODUCT INFORMATION



## MRT6160

Item No. 44274

**CAS Registry No.:** 3050683-38-7  
**Formal Name:** 3-[2-chloro-4'-(2-oxo-1(2H)-pyridinyl)[1,1'biphenyl]-3-yl]-2,6-piperidinedione  
**Synonym:** VAV1 Degradar-3  
**MF:** C<sub>22</sub>H<sub>17</sub>ClN<sub>2</sub>O<sub>3</sub>  
**FW:** 392.8  
**Purity:** ≥98%  
**Supplied as:** A 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

MRT6160 is supplied as a solid. A stock solution may be made by dissolving the MRT6160 in the solvent of choice, which should be purged with an inert gas. MRT6160 is slightly soluble (0.1-1 mg/ml) in DMSO.

### Description

MRT6160 is a degrader of proto-oncogene vav (VAV1).<sup>1</sup> It induces degradation of VAV1 with a half-maximal degradation concentration (DC<sub>50</sub>) value of 7 nM.

### Reference

1. Mcallister, L.A., Liardo, E., Ritzen, A., *et al.* Targeted degradation of VAV1. *Monte Rosa Therapeutics, Inc. WO2024/151547A1* (2024).

#### 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, 09/18/2025

#### 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