

# PRODUCT INFORMATION



## RAF265

Item No. 16991

**CAS Registry No.:** 927880-90-8  
**Formal Name:** 1-methyl-5-[[2-[5-(trifluoromethyl)-1H-imidazol-2-yl]-4-pyridinyl]oxy]-N-[4-(trifluoromethyl)phenyl]-1H-benzimidazol-2-amine

**Synonym:** CHIR265

**MF:** C<sub>24</sub>H<sub>16</sub>F<sub>6</sub>N<sub>6</sub>O

**FW:** 518.4

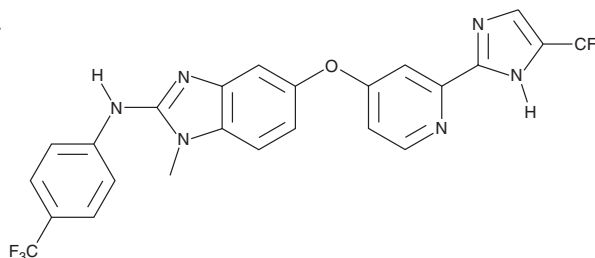
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 210, 266, 314 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

RAF265 is supplied as a crystalline solid. A stock solution may be made by dissolving the RAF265 in the solvent of choice, which should be purged with an inert gas. RAF265 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of RAF265 in these solvents is approximately 5, 10, and 2 mg/ml, respectively.

### Description

RAF265 is a dual inhibitor of mutant B-Raf<sup>V600E</sup> and vascular endothelial growth factor receptor 2 (EC<sub>50</sub>s = 0.14 and 0.19 μM, respectively, in cell-based assays).<sup>1</sup> It can also inhibit wild-type B-Raf, platelet-derived growth factor receptor, colony-stimulating factor 1R, RET, c-KIT, SRC, and STE20 with IC<sub>50</sub> values ranging from 20-100 nM.<sup>2</sup> RAF265 can inhibit proliferation of melanoma and colorectal cancer cell lines with active BRAF mutations as well as suppress tumor growth in melanoma xenograft models bearing B-Raf<sup>V600E</sup> mutations.<sup>2</sup>

### References

- Huang, T., Karsy, M., Zhuge, J., *et al.* B-Raf and the inhibitors: From bench to bedside. *J. Hematol. Oncol.* **6**, 1-9 (2013).
- Su, Y., Vilgelm, A.E., Kelley, M.C., *et al.* RAF265 inhibits the growth of advanced human melanoma tumors. *Clin. Cancer Res.* **18(8)**, 2184-2198 (2012).

#### 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, 11/04/2022

#### 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