# **PRODUCT** INFORMATION



## Vapendavir

Item No. 30546

CAS Registry No.:	439085-51-5
Formal Name:	3-ethoxy-6-[2-[1-(6-methyl-3-pyridazinyl)-
	4-piperidinyl]ethoxy]-1,2-benzisoxazole
Synonym:	BTA798
MF:	$C_{21}H_{26}N_4O_3$
FW:	382.5 N
Purity:	≥95%
UV/Vis.:	λ <sub>max</sub> : 219, 251 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

Vapendavir is supplied as a crystalline solid. A stock solution may be made by dissolving the vapendavir in the solvent of choice, which should be purged with an inert gas. Vapendavir is soluble in organic solvents such as chloroform and dimethyl formamide. The solubility of vapendavir in these solvents is approximately 10 and 1 mg/ml, respectively.

#### Description

Vapendavir is a capsid-binding antiviral agent.<sup>1,2</sup> It binds to a pocket in the human rhinovirus (HRV) strain HRV2 VP1 capsid protein and inhibits the cytopathic effects of HRV2 and HRV14 in HeLa Ohio cells ( $EC_{50}s = 1$  and 5 ng/ml, respectively), as well as in a panel of 39 HRV clinical isolates (median EC<sub>50</sub> = 7.3 ng/ml).<sup>1</sup> Vapendavir inhibits replication in a panel of 21 enterovirus 71 (EV71) isolates with an average  $EC_{50}$  value of 0.7  $\mu$ M.<sup>2</sup>

#### References

- 1. Feil, S.C., Hamilton, S., Krippner, G.Y., et al. An orally available 3-ethoxybenzisoxazole capsid binder with clinical activity against human rhinovirus. ACS Med. Chem. Lett. 3(4), 303-307 (2012).
- Tijsma, A., Franco, D., Tucker, S., et al. The capsid binder vapendavir and the novel protease inhibitor SG85 2. inhibit enterovirus 71 replication. Antimicrob. Agents Chemother. 58(11), 6990-6992 (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.

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, 11/16/2022

## 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