

PRODUCT INFORMATION



Tipranavir

Item No. 21028

CAS Registry No.: 174484-41-4

Formal Name: N-[3-[(1R)-1-[(6R)-5,6-dihydro-4-hydroxy-2-oxo-6-(2-phenylethyl)-6-propyl-2H-pyran-3-yl]propyl]phenyl]-5-(trifluoromethyl)-2-pyridinesulfonamide

Synonym: PNU-140690

MF: $C_{31}H_{33}F_3N_2O_5S$

FW: 602.7

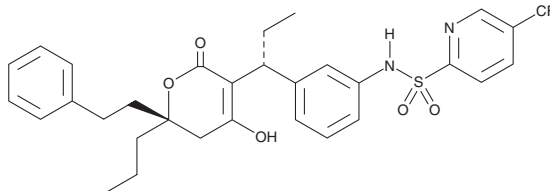
Purity: $\geq 95\%$

UV/Vis.: λ_{max} : 254 nm

Supplied as: A crystalline solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Tipranavir is supplied as a crystalline solid. A stock solution may be made by dissolving the tipranavir in the solvent of choice, which should be purged with an inert gas. Tipranavir is slightly soluble in chloroform and methanol.

Description

Tipranavir is a potent, orally bioavailable nonpeptidic HIV aspartyl protease inhibitor ($K_i = < 1$ nM) that prevents cleavage and thus activation of functional HIV cleaved proteins.^{1,2} Formulations containing tipranavir, in combination with other HIV treatments, are effective against multidrug resistant viruses and are used in the treatment of HIV infection progression in patients who have not responded well to initial treatments.³

References

1. Turner, S.R., Strohbach, J.W., Tommasi, R.A., et al. Tipranavir (PNU-140690): A potent, orally bioavailable nonpeptidic HIV protease inhibitor of the 5,6-dihydro-4-hydroxy-2-pyrone sulfonamide class. *J. Med. Chem.* **41**(18), 3467-3476 (1998).
2. Thaisrivongs, S., and Strohbach, J.W. Structure-based discovery of Tipranavir disodium (PNU-140690E): A potent, orally bioavailable, nonpeptidic HIV protease inhibitor. *Biopoly.* **51**(1), 51-58 (1999).
3. Llibre, J.M., Schapiro, J.M., and Clotet, B. Clinical implications of genotypic resistance to the newer antiretroviral drugs in HIV-1-infected patients with virological failure. *Clin. Infect. Dis.* **50**(6), 872-881 (2010).

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, 11/16/2022

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