

PRODUCT INFORMATION



THK5351

Item No. 45302

CAS Registry No.: 1707147-26-9

Formal Name: 1-fluoro-3-[[2-[6-(methylamino)-3-pyridinyl]-6-quinolinyl]oxy]-2S-propanol

MF: C₁₈H₁₈¹⁸FN₃O₂

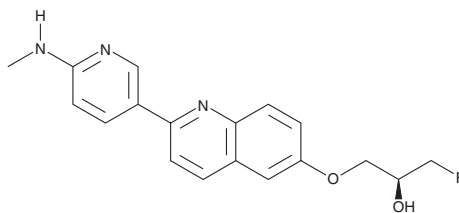
FW: 327.4

Purity: ≥95%

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

THK5351 is supplied as a solid. A stock solution may be made by dissolving the THK5351 in the solvent of choice, which should be purged with an inert gas. THK5351 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

THK5351 is a non-radiolabeled form of ¹⁸F-THK5351, which has been used as a PET tracer for selectively imaging tau in animal models of tauopathies.^{1,2}

References

1. Moreno-Gonzalez, I., Edwards, G.A., 3rd, Hasan, O., *et al.* Longitudinal assessment of tau-associated pathology by ¹⁸F-THK5351 pet imaging: A histological, biochemical, and behavioral study. *Diagnostics (Basel)* **11**(10), 1874 (2016).
2. Harada, R., Okamura, N., Furumoto, S., *et al.* ¹⁸F-THK5351: A Novel pet radiotracer for imaging neurofibrillary pathology in Alzheimer disease. *J. Nucl. Med.* **57**(2), 208-214 (2016).

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, 04/14/2026

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