

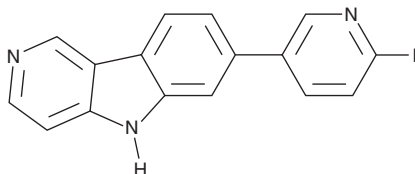
PRODUCT INFORMATION



T807

Item No. 42289

CAS Registry No.: 1415379-56-4
Formal Name: 7-(6-fluoro-3-pyridinyl)-5H-pyrido[4,3-b]indole
MF: $C_{16}H_{10}FN_3$
FW: 263.3
Purity: $\geq 98\%$
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

T807 is supplied as a solid. A stock solution may be made by dissolving the T807 in the solvent of choice, which should be purged with an inert gas. T807 is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

T807 is a tau-binding agent.¹ It selectively binds to recombinant tau ($K_d = 0.2$ and 16.8 nM for high- and low-affinity tau binding sites) over 83 receptors, transporters, enzymes, and ion channels in a panel at $10\text{ }\mu\text{M}$ but does not bind to monoamine oxidase A (MAO-A) and MAO-B (K_d s = 1.6 and 21 nM, respectively). T807 binds to native tau fibrils from patients with Alzheimer's disease or progressive supranuclear palsy (PSP; K_d s = 0.3 nM for both for high-affinity tau binding sites; K_d s = 30 and 70 nM, respectively, for low-affinity tau binding sites) but also binds to brain homogenates from patients without a tau-related disease (K_d s = 0.6 and 52.5 nM for high- and low-affinity tau binding sites, respectively) due to its off-target binding to MAO-A and MAO-B. A radiolabeled form of T807, ^{18}F -T807, has been used in the imaging of tau in animal models of tau pathology *in vivo* and in postmortem human brain tissue from patients with Alzheimer's disease.²

References

1. Vermeiren, C., Motte, P., Viot, D., *et al.* The tau positron-emission tomography tracer AV-1451 binds with similar affinities to tau fibrils and monoamine oxidases. *Mov. Disord.* **33**(2), 273-281 (2018).
2. Brendel, M., Yousefi, B.H., Blume, T., *et al.* Comparison of ^{18}F -T807 and ^{18}F -THK5117 PET in a mouse model of tau pathology. *Front. Aging Neurosci.* **10**, 174 (2018).

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, 11/22/2024

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