

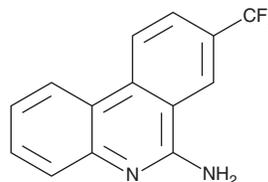
# PRODUCT INFORMATION



## 6-Amino-8-trifluoromethylphenanthridine

Item No. 26419

**CAS Registry No.:** 651055-83-3  
**Formal Name:** 8-(trifluoromethyl)-6-phenanthridinamine  
**Synonym:** 6A-8tFP  
**MF:** C<sub>14</sub>H<sub>9</sub>F<sub>3</sub>N<sub>2</sub>  
**FW:** 262.2  
**Purity:** ≥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

6-Amino-8-trifluoromethylphenanthridine (6A-8tFP) is supplied as a solid. A stock solution may be made by dissolving the 6A-8tFP in the solvent of choice, which should be purged with an inert gas. 6A-8tFP is soluble in organic solvents such as ethanol and DMSO.

### Description

6A-8tFP is an antiprion agent and a derivative of 6-aminophenanthridine (Item No. 26420).<sup>1</sup> It inhibits protein folding activity of the ribosome (PFAR) when used at a concentration of 150 μM.<sup>2</sup> 6A-8tFP directly competes with protein substrates for the ribosomal active site.

### References

1. Bach, S., Talarek, N., Andrieu, T., *et al.* Isolation of drugs active against mammalian prions using a yeast-based screening assay. *Nat. Biotechnol.* **21(9)**, 1075-1081 (2003).
2. Banerjee, D., Vovusha, H., Pang, Y., *et al.* Spectroscopic and DFT studies on 6-Aminophenanthridine and its derivatives provide insights in their activity towards ribosomal RNA. *Biochimie* **97**, 194-199 (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.

#### 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, 12/20/2022

#### 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