

TTFA

Mitochondrial Complex II inhibitor

TTFA blocks the respiratory chain complex II causing inhibition of mitochondrial respiration. Respiratory chain complex II inhibition is caused via binding of TTFA to two ubiquinone binding sites, Qp and Qd. Inhibition of Complex II by TFA has been shown to cause a delay in overall cell cycle progression leading to oxidative stress. TTFA also was found to inhibit porcine liver carboxylesterase.

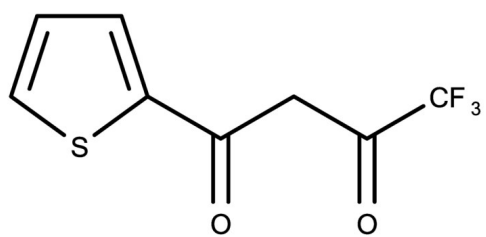
Ordering Information

[Order Online »](#)

ENZ-CHM497-0500	500mg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Thenoyltrifluoroacetone, 4,4,4-Trifluoro-1-(2-thienyl)-1,3-butanedione

Appearance Pale yellow solid.

CAS 326-91-0

Couple Type Inhibitor

Formula $C_8H_5F_3O_2S$

Identity Determined by NMR.

MW 222.18

Purity ≥98% (TLC)

Solubility Soluble in DMSO (25 mg/ml) or ethanol (25 mg/ml).

Last modified: February 27, 2025



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)