

TOFA

Acetyl-CoA inhibitor

Hypolipidemic agent. Potent, reversible and competitive inhibitor of acetyl-coenzyme A carboxylase (Ac-CoA carboxylase; ACC), a key enzyme involved in fatty acid biosynthesis. It blocks apoptosis induced by the fatty acid synthase inhibitor C75 (Prod. No. ALX-270-286) and stimulates neurite outgrowth and neuronal differentiation in rat pheochromocytoma cells.

Citations: 11

[View Online »](#)

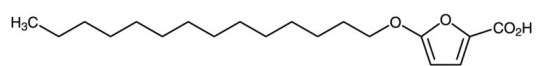
Ordering Information

[Order Online »](#)

| | |
|----------------|------|
| BML-EI321-0010 | 10mg |
| BML-EI321-0050 | 50mg |

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

| | |
|-------------------|---|
| Use/Stability | As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C. |
| Long Term Storage | -20°C |
| Shipping | Blue Ice |

Regulatory Status

RUO - Research Use Only

Product Details

| | |
|------------------|--|
| Alternative Name | 5-(Tetradecyloxy)-2-furoic acid, RMI-14514 |
| Appearance | Off-white solid. |
| CAS | 54857-86-2 |
| Couple Target | Acetyl-CoA |
| Couple Type | Inhibitor |
| Formula | $C_{19}H_{32}O_4$ |
| MW | 324.5 |
| MeltingPoint | 108-110°C |
| Purity | ≥99% (TLC) |
| Solubility | Soluble in dimethyl formamide, DMSO (10mg/ml), 100% ethanol or methanol. Sparingly soluble in aqueous buffers. |



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-
fr@enzolifesciences.com](mailto:info-fr@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)