

Sobetirome

thyroid hormone receptor TR β 1 agonist

Sobetirome, also known as GC-1, is a selective thyromimetic drug that mimics the action of thyroid hormones. It binds preferentially to the thyroid hormone receptor TR β 1 over TR α 1, which helps to minimize the typical side effects associated with thyroid hormone therapies.

Ordering Information

[Order Online »](#)

ENZ-CHM329-0005	5mg
ENZ-CHM329-0025	25mg

Manuals, SDS & CofA

[View Online »](#)

- GMP format available

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.Powder: Stable at -20°C for 3 years or +4°C for 2 years; In solvent: Stable at -80°C for 2 years or -20°C for 1 year.
Handling	Before opening the vial, please centrifuge to gather the compound at the bottom of the vial.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	GC-1, QRX-43, 2-[4-[[4-hydroxy-3-(1-methylethyl)phenyl]methyl]-3,5-dimethylphenoxy]-acetic acid
Appearance	White to light yellow solid.
CAS	211110-63-3
Couple Target	Thyroid hormone receptor
Couple Type	Ligand
Formula	C ₂₀ H ₂₄ O ₄
Identity	Identity determined by LCMS.
MW	328.4
MeltingPoint	122-124°C
Purity	≥99% (LCMS)
Solubility	Soluble in DMSO (100 mg/ml).



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)