

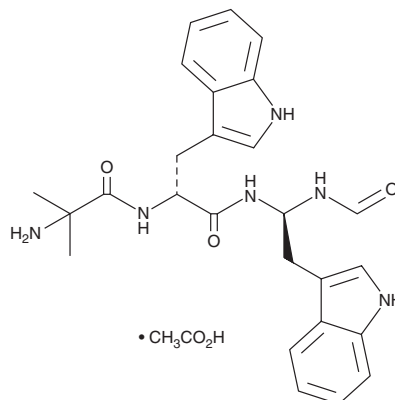
PRODUCT INFORMATION



Macimorelin (acetate)

Item No. 41886

CAS Registry No.: 945212-59-9
Formal Name: 2-methylalanyl-N-[(1R)-1-(formylamino)-2-(1H-indol-3-yl)ethyl]-D-typtophanamide, acetate
Synonyms: AEZS-130, Aib-D-Trp-D-gTrp-CHO, Aib-D-Trp-D-gTrp-formyl, D-87575, EP 1572, JMV1843
MF: $C_{26}H_{30}N_6O_3 \cdot C_2H_4O_2$
FW: 534.6
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

Macimorelin (acetate) is supplied as a solid. A stock solution may be made by dissolving the macimorelin (acetate) in the solvent of choice, which should be purged with an inert gas. Macimorelin (acetate) is soluble (≥ 10 mg/ml) in ethanol and DMSO.

Description

Macimorelin is a growth hormone secretagogue and agonist of growth hormone secretagogue receptor 1a (GHS-R1a).¹ It binds to GHS-R1a ($IC_{50} = 123$ nM for the human receptor). Oral administration of macimorelin (1 mg/kg) increases plasma growth hormone levels in dogs. It decreases gastric acid secretion in pylorus-ligated rats when administered intracerebroventricularly at doses of 0.01 or 1 nmol/animal.² Macimorelin (10 mg/kg) increases acute food intake and body weight in mice.³ It decreases the number and total duration of seizures in mice induced with dopamine D_1 receptor kindling using SKF 81297.⁴ Formulations containing macimorelin have been used in the diagnosis of adult growth hormone deficiency.

References

1. Guerlavais, V., Boeglin, D., Mousseaux, D., *et al.* New active series of growth hormone secretagogues. *J. Med. Chem.* **46**(7), 1191-1203 (2003).
2. Sibilia, V., Muccioli, G., Deghenghi, R., *et al.* Evidence for a role of the GHS-R_{1a} receptors in ghrelin inhibition of gastric acid secretion in the rat. *J. Neuroendocrinol.* **18**(2), 122-128 (2006).
3. Holubová, M., Spolcová, A., Demianová, Z., *et al.* Ghrelin agonist JMV 1843 increases food intake, body weight and expression of orexigenic neuropeptides in mice. *Physiol. Res.* **62**(4), 435-444 (2013).
4. Buckinx, A., Van Den Herrewegen, Y., Pierre, A., *et al.* Differential effects of a full and biased ghrelin receptor agonist in a mouse kindling model. *Int. J. Mol. Sci.* **20**(10), 2480 (2019).

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, 10/16/2024

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