

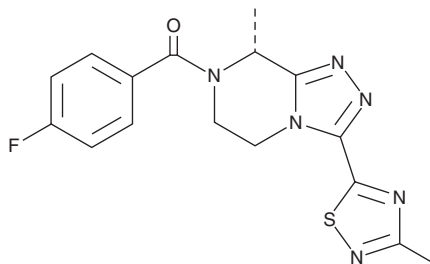
PRODUCT INFORMATION



Fezolinetant

Item No. 39523

CAS Registry No.: 1629229-37-3
Formal Name: [(8R)-5,6-dihydro-8-methyl-3-(3-methyl-1,2,4-thiadiazol-5-yl)-1,2,4-triazolo[4,3-a]pyrazin-7(8H-yl)](4-fluorophenyl)-methanone
Synonyms: AS 3472693-00, ESN364
MF: C₁₆H₁₅FN₆OS
FW: 358.4
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

Fezolinetant is supplied as a solid. A stock solution may be made by dissolving the fezolinetant in the solvent of choice, which should be purged with an inert gas. Fezolinetant is soluble in the organic solvent DMSO.

Description

Fezolinetant is a neurokinin-3 (NK₃) receptor antagonist (IC₅₀ = 20 nM).¹ It decreases plasma levels of luteinizing hormone (LH) in castrated cynomolgus monkeys when administered at a dose of 5 mg/kg. Fezolinetant (10 mg/kg) prevents increases in body weight and food intake but does not affect uterine weight or plasma levels of estradiol (17β-estradiol; Item No. 10006315), follicle-stimulating hormone (FSH), progesterone (Item No. 15876), or testosterone in ovariectomized rats.² It inhibits increases in skin temperature but not core temperature, indicating regulation of hot flash-like symptoms, in ovariectomized rats when administered at doses of 3 or 10 mg/kg. Formulations containing fezolinetant have been used in the treatment of moderate to severe vasomotor symptoms due to menopause.

References

1. Hoveyda, H.R., Fraser, G.L., Dutheuil, G., *et al.* Optimization of novel antagonists to the neurokinin-3 receptor for the treatment of sex-hormone disorders (Part II). *ACS Med. Chem. Lett.* **6**(7), 736-740 (2015).
2. Tahara, A., Takamatsu, H., Ohtake, A., *et al.* Effects of neurokinin 3 receptor antagonist fezolinetant on hot flash-like symptoms in ovariectomized rats. *Eur. J. Pharmacol.* **905**, 174207 (2021).

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, 10/16/2023

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