

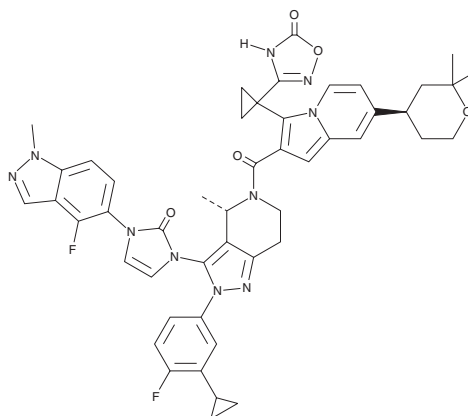
PRODUCT INFORMATION



ECC5004

Item No. 41324

CAS Registry No.: 3011682-51-9
Formal Name: 3-[1-[2-[[[4S]-2-(3-cyclopropyl-4-fluorophenyl)-3-(4-fluoro-1-methyl-1H-indazol-5-yl)-2,3-dihydro-2-oxo-1H-imidazol-1-yl]-2,4,6,7-tetrahydro-4-methyl-5H-pyrazolo[4,3-c]pyridin-5-yl] carbonyl]-7-[(4R)-tetrahydro-2,2-dimethyl-2H-pyran-4-yl]-3-indoliziny] cyclopropyl]-1,2,4-oxadiazol-5(2H)-one



Synonym: AZD 5004
MF: C₄₈H₄₆F₂N₁₀O₅
FW: 880.9
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

ECC5004 is supplied as a solid. A stock solution may be made by dissolving the ECC5004 in the solvent of choice, which should be purged with an inert gas. ECC5004 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

ECC5004 is an agonist of glucagon-like peptide 1 receptor (GLP-1R).¹ It induces cAMP accumulation in HEK293 cells expressing human GLP-1R with an EC₅₀ value less than 20 nM.

References

1. Ren, Z., Sun, X., Zhou, J., *et al.* Substituted imidazoles as GLP-1 receptor agonists. *Eccogene (Shanghai) Co., Ltd.* US 11,584,751 B1 (2023).

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, 10/15/2024

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