PRODUCT INFORMATION



Herboxidiene

Item No. 25136

| CAS Registry No.: | 142861-00-5 |
|------------------------------------------------------------------------------------------------------------------------------------|------------------------------------------------|
| Formal Name: | 5,6-anhydro-6-C-[(2S,3E,5E)-6- |
| | [(2S,3S,6R)-6-(carboxymethyl)tetrahydro- |
| | 3-methyl-2H-pyran-2-yl]-2-methyl-3,5- |
| | heptadien-1-yl]-1,4,7-trideoxy-4-methyl- |
| | 3-O-methyl-L-glycero-L-gluco-heptitol |
| Synonyms: | GEX1A, Tan 1609 он |
| MF: | C ₂₅ H ₄₂ O ₆ |
| FW: | 438.6 |
| Purity: | ≥70% |
| Supplied as: | A solid Ö |
| Storage: | -20°C |
| Stability: | ≥4 years |
| Item Origin: | Bacterium/Streptomyces sp. |
| Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis. | |

Laboratory Procedures

Herboxidiene is supplied as a solid. A stock solution may be made by dissolving the herboxidiene in the solvent of choice. Herboxidiene is soluble in organic solvents such as ethanol, methanol, DMSO, and dichloromethane, which should be purged with an inert gas.

Description

Herboxidiene is a polyketide originally isolated from S. chromofuscus that has diverse biological activities.¹⁻⁵ It inhibits growth of HeLa S3, SK-MEL-2, PC3, A549, and EBC-1 cells with GI₅₀ values ranging from 7.4 to 62 nM.1 Herboxidiene is cytostatic against human umbilical vein endothelial cells (HUVECs; $(IC_{50} = 26 \text{ nM}))$ and inhibits VEGF-induced invasion and tube formation of serum-starved HUVECs in a concentration-dependent manner, indicating antiangiogenic activity.² Herboxidiene (0.05 μ M) inhibits HIF-1a mRNA splicing and reduces HIF-1a protein levels in HepG2 cells grown under hypoxic conditions. It also inhibits splicing of $p27^{Kip}$ mRNA in HeLa cells in a concentration-dependent manner via interaction with the SAP155 subunit of the SF3b complex.³ Herboxidiene (0.1 and 1 μ M) increases LDL receptor promoter-driven transcription in a cell-based reporter assay.⁴

References

- 1. Imaizumi, T., Nakagawa, H., Hori, R., et al. J. Antibiot. (Tokyo) 70(5), 675-679 (2017).
- 2. Jung, H.J., Kim, Y., Shin, J.Y., et al. Arch. Pharm. Res. 38(9), 1728-1735 (2015).
- 3. Hasegawa, M., Miura, T., Kuzuya, K., et al. ACS Chem Biol. 6(3), 229-233 (2011).
- 4. Koguchi, Y., Nishio, M., Kotera, J., et al. J. Antibiot. (Tokyo) 50(11), 970-971 (1997).
- 5. Miller-Wideman, M., Makkar, N., Tran, M.G.B., et al. J. Antibiot. (Tokyo) 45(6), 914-921 (1992).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFFTY 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

uyer 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, 08/28/2023

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