

# PRODUCT INFORMATION



MA-0204

Item No. 26025

**CAS Registry No.:** 2095128-17-7  
**Formal Name:** 3R-methyl-6-[2-[[5-methyl-2-[4-(trifluoromethoxy)phenyl]-1H-imidazol-1-yl]methyl]phenoxy]-hexanoic acid

**MF:** C<sub>25</sub>H<sub>27</sub>F<sub>3</sub>N<sub>2</sub>O<sub>4</sub>

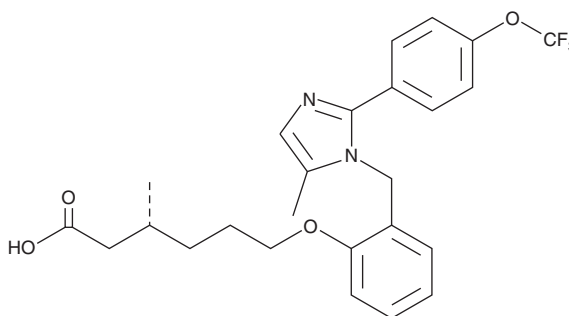
**FW:** 476.5

**Purity:** ≥95%

**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

MA-0204 is supplied as a solid. A stock solution may be made by dissolving the MA-0204 in the solvent of choice, which should be purged with an inert gas. MA-0204 is soluble in methanol and DMSO.

## Description

MA-0204 is an agonist of peroxisome proliferator-activated receptor  $\delta$  (PPAR $\delta$ ).<sup>1</sup> It selectively induces reporter gene expression in CV-1 cells expressing human PPAR $\delta$  (EC<sub>50</sub> = 0.4 nM) over CV-1 cells expressing human PPAR $\alpha$  or PPAR $\gamma$  (EC<sub>50</sub>s = 6,990 and >10,000 nM, respectively). MA-0204 increases the expression of mRNA encoding angiopoietin-related protein 4 (ANGPTL4), carnitine palmitoyltransferase 1 (CPT1), and pyruvate dehydrogenase kinase 4 (PDHK4) in primary myoblasts from a patient with Duchenne muscular dystrophy (DMD) in a concentration-dependent manner. *In vivo*, MA-0204 (30 or 100 mg/kg) increases the expression of mRNA encoding Pdhk4 in the skeletal muscle of mice.

## Reference

- Scholtens, P.A.M.J., Oozeer, R., Martin, R., *et al.* The early settlers: Intestinal microbiology in early life. *Annu. Rev. Food Sci. Technol.* **3**, 425-447 (2012).

### 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, 02/22/2024

## 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