

Discovery Life Sciences
6 Henshaw Street
Woburn, MA 01801
Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

4'-Hydroxydiclofenac-[13C6]

Catalog Number: 451006

Storage Conditions: STORE AT -20°C

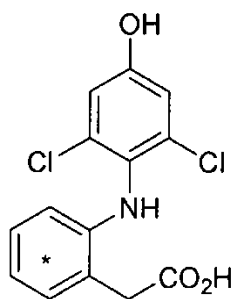
Lot Number: 2406250

Date Released:.....2024 June

Expiration Date:.....2026 June

Amount: 1 mg

Structure:



Molecular Formula: $^{13}\text{C}_6^{12}\text{C}_8\text{H}_{11}\text{Cl}_2\text{NO}_3$

Formula Weight: 318.15

CAS Number: 64118-84-9 (unlabeled)

Purity: $\geq 99\%$ by HPLC chromatography with absorbance detection.

Identity:

HPLC Chromatography: Conforms to standard

Absorbance Spectrum: Conforms to standard

$^1\text{H-NMR}$: Conforms to structure

Mass Spectrum: Conforms to structure

Use: Dissolve in ACN or MeOH at a concentration of 1 mM.

4'-Hydroxydiclofenac-[13C6]

4'-Hydroxydiclofenac-[13C6] is the reference metabolite for the Diclofenac-4-Hydroxylase Assay for CYP2C9. This reagent can be used as an internal standard for enzymatic reactions that are analyzed by LCMS.

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© DLS

Discovery Life Sciences
6 Henshaw Street
Woburn, MA 01801
Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

SAFETY RECOMMENDATIONS:

Safety assessment indicates this product is hazardous; refer to SDS [Safety Data Sheet] for safety information.

Handle in accordance with good industrial hygiene and laboratory safety practices.



01 July 2024

Quality Assurance

Date

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© DLS