

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

## 4'-Hydroxydiclofenac-[13C6]

**Catalog Number:** 451006

**Storage Conditions:** STORE AT -20°C

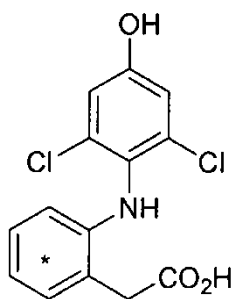
**Lot Number:** 2408297

**Date Released:** 2024 September

**Expiration Date:**.....2025 January

**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:** 99 % by HPLC chromatography with absorbance detection .

**Identity:**

*HPLC Chromatography:* Conforms to standard

*Absorbance Spectrum:* Conforms to standard

*<sup>1</sup>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](mailto: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.



---

**Quality Assurance**

10 September 2024

---

**Date**

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

© DLS