

---

## PARP Inhibitor XIV

(Catalog No. M77000)

---

### Alternative Names

4-(1-Methyl-1H-pyrrol-2-ylmethylene)-4H-isoquinoline-1,3-dione

### Description

A cell-permeable isoquinolinedione compound that potently inhibits PARP-1 activity in a NAD<sup>+</sup>-competitive and reversible manner. Affects PPAR-2 activity with lower potency. Selectively inhibits PARP-1-mediated cellular PAR formation and is of little effect against PAR formation in fibroblasts derived from *parp-1*(-/-) mice. Its in vivo application is limited

### CAS Number

1104546-89-5

### Chemical Formula

C<sub>15</sub>H<sub>12</sub>N<sub>2</sub>O<sub>2</sub>

### Molecular Weight

252.3

### Size

10 mg

### Purity

> 97%

### Solubility

Soluble in DMSO

### Storage Temp

-20°C

### Ordering Information:

Products	Size	Cat. No.
PARP Inhibitor XIV	10 mg	M77000 -1

*This product is for research purposes only. Not intended for use in diagnostic procedures.*