



**EAGLE**  
BIOSCIENCES

**Anti DMPO Rabbit Polyclonal Antiserum**

Catalog Number: DMP00-A100

Aliquot: 100  $\mu$ l

Storage: -70°C

**PRODUCT SPECIFICATION SHEET**

<b>Immunogen:</b>	5,5-dimethyl-1-pyrroline-N-oxide (DMPO) – octanoic acid adduct conjugated to ovalbumin.
<b>Concentration:</b>	unpurified antiserum
<b>Aliquot:</b>	100 $\mu$ L
<b>Applications:</b>	Western Blot- 1:5,000 ELISA- 1:5,000 to 1:10,000 IHC- 1:1,000 subject to optimization
<b>Storage and Stability:</b>	At least one year at -70°C. AVOID REPEATED FREEZE THAW CYCLES.

*Product Developed and Manufactured in the USA*

---

EAGLE BIOSCIENCES, INC.  
20A NW Blvd., Suite 112  
Nashua, NH 03063

Phone: 617-419-2019  
Fax: 617-419-1110  
Email: [info@EagleBio.com](mailto:info@EagleBio.com)  
[www.EagleBio.com](http://www.EagleBio.com)