



**EAGLE**  
BIOSCIENCES

**Anti-Human Estrogen Receptor  $\beta$   
(ER $\beta$ )**

Product Number: ERB00-A100C

Aliquot: 0.1 mg

Storage: -20° C

**PRODUCT SPECIFICATION SHEET**

<b>IMMUNOGLOBULIN:</b>	Polyclonal Chicken IgY
<b>IMMUNOGEN:</b>	Recombinant Human VEGF
<b>PREDICTED SIZE:</b>	23 kDa (monomer), 37 kDa (dimer)
<b>CONCENTRATION:</b>	1.0 mg/mL
<b>SPECIFICITY:</b>	Human VEGF, not tested on other species.
<b>STORAGE:</b>	Store at -20° C. Avoid repeating freeze/thaw cycles, as it may result in loss of activity. Store at 4° C for short term.
<b>PURITY:</b>	IgY
<b>WORKING DILUTION:</b>	Suggested concentration for Western blots is 1.0 $\mu$ g/mL. Not tested for IHC or ELISA.

*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)