



EAGLE
BIOSCIENCES

Anti-6-keto-PGF1 α

Product Number: KET00-A1000

Aliquot: 1 mL

Storage: -70° C

PRODUCT SPECIFICATION SHEET

Contents: Sheep polyclonal antiserum specific for 6-keto-PGF1 alpha in phosphate buffered saline with 0.1% sodium azide.

Quantity: 1 mL. Sufficient for at least 100 tubes in conventional liquid phase radioimmunoassay procedures.

Use: Thaw antiserum at room temperature. Make a 1:25 working dilution of the reconstituted antiserum in PBS plus 0.1% gelatin (i.e. assay buffer). This permits addition of 200 μ L per assay tube.

Sensitivity: The suggested dilution will result in >30% binding of tracer in a liquid phase RIA with 0.2 ml of first antibody added to a final volume of 1.0 mL.

Specificity:

<u>Hapten:</u>	<u>Reactivity (%)</u>
PGF2	2.10
6-ketoPGF1	100.00
thromboxane B2	0.07
13,14-dihydro-15-keto PG2	0.03
13,14-dihydro-6.15-diketo PGF2	0.02
PGE2	0.05
PGD2	0.03
PGE1	0.06
PGE2	0.05
6-ketoPGE1	8.60
PGF1	7.90
arachidonic acid	0.002

References:

J.L. Fahey, et al., Amer. Soc. Microbiol. pp. 685-691 (1986)
W.S. Powell, Methods in Enzymology **86**;467-476.

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
www.EagleBio.com