



**EAGLE**  
BIOSCIENCES

## Anti-Thromboxane B<sub>2</sub> Polyclonal Antibody

Product Number: TBB00-A1000A

Aliquot: 1.0 mL

Storage: -80°C

### PRODUCT SPECIFICATION SHEET

<b>DESCRIPTION:</b>	Rabbit antisera specific for thromboxane B <sub>2</sub> .												
<b>CONCENTRATION:</b>	Not Determined												
<b>STORAGE:</b>	-80°C. <b>AVOID REPEATED FREEZE THAWS.</b>												
<b>STORAGE BUFFER:</b>	PBS with 0.1% sodium azide												
<b>IMMUNOGEN:</b>	Thromboxane B <sub>2</sub>												
<b>PURIFICATION:</b>	Not Purified												
<b>APPLICATIONS:</b>	Suitable for use in RIA. May also have utility in ELISA.												
<b>SENSITIVITY:</b>	50 pg of unlabeled TXB <sub>2</sub> will result in 50% inhibition of radiolabeled ligand binding in an RIA system.												
<b>SPECIES SPECIFICITY:</b>	Not species specific. May not exhibit utility when using in a rabbit experimental model.												
<b>REFERENCE:</b>	Hamberg M, Svensson J and Samuelsson B (1975). <i>Proc. Natl. Acad. Sci. USA</i> <u>72</u> :2994-2998. Oates JA, FitzGerald GA, Branch A, Jackson EK, Knapp HR and Roberts LJ (1988). <i>N. Engl. J. Med.</i> <u>319</u> :689-698. Powell WS (1982). <i>Meth. Enzymol.</i> <u>86</u> :467-476.												
<b>CROSS-REACTIVITY:</b>	<table><thead><tr><th>HAPTEN</th><th>REACTIVITY</th></tr></thead><tbody><tr><td>TXB<sub>2</sub></td><td>100%</td></tr><tr><td>PGD<sub>2</sub></td><td>&lt;1%</td></tr><tr><td>PGE<sub>2</sub></td><td>&lt;1%</td></tr><tr><td>PGF<sub>2α</sub></td><td>&lt;1%</td></tr><tr><td>6,keto-PGF<sub>1α</sub></td><td>&lt;1%</td></tr></tbody></table>	HAPTEN	REACTIVITY	TXB <sub>2</sub>	100%	PGD <sub>2</sub>	<1%	PGE <sub>2</sub>	<1%	PGF <sub>2α</sub>	<1%	6,keto-PGF <sub>1α</sub>	<1%
HAPTEN	REACTIVITY												
TXB <sub>2</sub>	100%												
PGD <sub>2</sub>	<1%												
PGE <sub>2</sub>	<1%												
PGF <sub>2α</sub>	<1%												
6,keto-PGF <sub>1α</sub>	<1%												

*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)