



**EAGLE**  
BIOSCIENCES

## Anti-Protein Kinase C $\beta$ II

Catalog Number: PK200-A1000

Aliquot: 1000  $\mu$ L

Storage: -70°C

### PRODUCT SPECIFICATION SHEET

<b>DESCRIPTION:</b>	This polyclonal antibody is produced in rabbit and is provided as antiserum specific for Protein Kinase C $\beta$ II (PKC $\beta$ II).
<b>MOLECULAR WEIGHT:</b>	77 kDa Calculated
<b>CONCENTRATION:</b>	Not Determined
<b>ISOTYPE:</b>	NA
<b>STORAGE BUFFER:</b>	Phosphate Buffered Saline; 0.05% Sodium Azide
<b>STORAGE:</b>	-70°C. AVOID REPEATED FREEZE THAW CYCLES. Has demonstrated stability at 4 °C for greater than 1 month.
<b>IMMUNOGEN:</b>	KLH conjugated synthetic peptide corresponding to amino acids 661- 673 of PKC $\beta$ II. <b>Sequence:</b> <sup>661</sup> F-V-N-S-E-F-L-K-P-E-V-K-S <sup>673</sup>
<b>PURIFICATION:</b>	NA
<b>APPLICATIONS:</b>	Western Blotting: 1:100 Dilution
<b>SPECIES SPECIFICITY:</b>	Human, mouse and rat among other mammalian species.
<b>REFERENCE:</b>	JBC (1993) <b>268</b> :20110-20115

*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)