

noassay. It can be used as a ubiquitous antiserum for the determination of plasma or in a range of fish species.

<b>References:</b>	Degger B <i>et al</i> (2000) <i>General and Comparative Endocrinology</i> <b>117</b> , 340-343
<b>Antigen:</b>	Barramundi ( <i>Lates calcarifer</i> ) IGF-I
<b>Species:</b>	Rabbit antiserum
<b>Specificity:</b>	Cross-reacts 100% with Barramundi IGF-I, Bream IGF-I ( <i>Pseudorasbora parva</i> ), Salmon and Trout IGF-I ( <i>Onchorynchus</i> and <i>Salmo</i> spp), Tilapia IGF-I ( <i>Oreochromis mossambicus</i> ) and Tuna IGF-I ( <i>Thunnus maccoyii</i> ). Cross-reactivity to human IGF-I and IGF-II is < 1%. No cross reactivity to salmon insulin.
<b>Form:</b>	Lyophilized powder
<b>Storage/ Stability:</b>	At least 2 years at 2 - 4°C (lyophilized). After reconstitution store at -20°C or -80°C Avoid freeze-thaw cycles
<b>Reconstitution:</b>	Dissolve the vial contents in 500µl RIA buffer (30mM sodium phosphate pH 7.5 containing 10mM EDTA, 0.2g/l potassium sulphate, 0.5ml Tween 20, 0.25g/l sodium azide) to give an antiserum stock of 1:50.
<b>Application and Titre:</b>	Each application and titre should be determined in house but as a guideline - Determination of IGF-I or IGF-II in a Range of Species by Radioimmunoassay (RIA). At a dilution of 1:30,000 the antiserum binds 20% of freshly prepared [ <sup>125</sup> I]- barramundi IGF-I (~80Ci/g). Half-maximal inhibition (ED <sub>50</sub> ) of binding is produced by 1.7 ng/ml of barramundi IGF-I.

## Product Codes and Pricing

**Barramundi IGF-I  
Antiserum**

**10 µl**

**PAF1**

**Related Products:**

Barramundi IGF-I

Bream IGF-I

Salmon/Trout IGF-I

Tuna IGF-I