

# Pierce® Avidin

21121 21128

0420.3

Number	Description
21121	Pierce Avidin (salt free), 10mg
21128	Pierce Avidin (salt free), 20mg

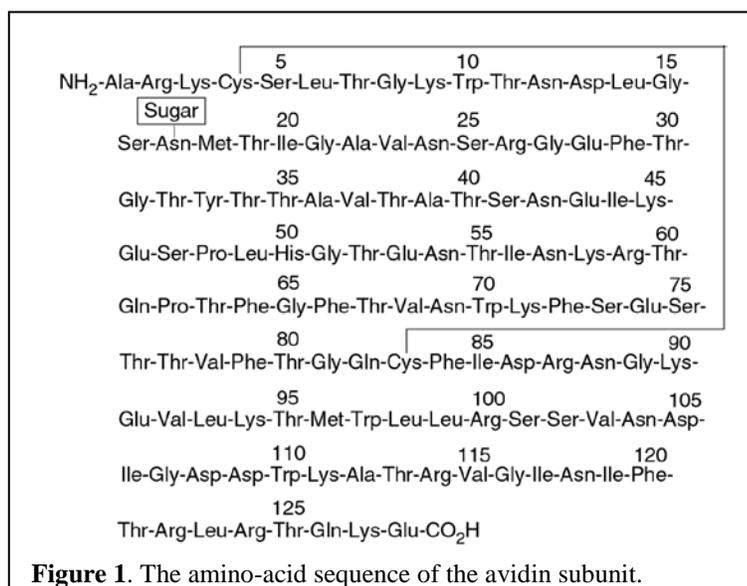
Specific Activity: 11-15µg biotin bound/mg protein

**Storage:** Upon receipt store product at -20°C. Product is shipped at ambient temperature.

## Introduction

Avidin is a basic glycoprotein present in the egg white and tissues of birds, reptiles and amphibia. This protein contains four identical subunits consisting of 128 amino acids (Figure 1) and has a combined molecular mass of 67-68K daltons.<sup>1</sup> Avidin has an isoelectric point of pH 10-10.5, is soluble in aqueous solutions and is stable over wide pH and temperature ranges.<sup>2</sup>

Thermo Scientific Pierce Avidin is a purified avidin protein. The avidin/biotin interaction is the strongest known noncovalent, biological interaction ( $K_a=10^{15} \text{ M}^{-1}$ ) between protein and ligand. The avidin/biotin complex is not significantly affected by pH values between 2 and 13 or by guanidine-HCl concentrations up to 8M at neutral pH.



This product ("Product") is warranted to operate or perform substantially in conformance with published Product specifications in effect at the time of sale, as set forth in the Product documentation, specifications and/or accompanying package inserts ("Documentation") and to be free from defects in material and workmanship. Unless otherwise expressly authorized in writing, Products are supplied for research use only. No claim of suitability for use in applications regulated by FDA is made. The warranty provided herein is valid only when used by properly trained individuals. Unless otherwise stated in the Documentation, this warranty is limited to one year from date of shipment when the Product is subjected to normal, proper and intended usage. This warranty does not extend to anyone other than the original purchaser of the Product ("Buyer").

**No other warranties, express or implied, are granted, including without limitation, implied warranties of merchantability, fitness for any particular purpose, or non infringement. Buyer's exclusive remedy for non-conforming Products during the warranty period is limited to replacement of or refund for the non-conforming Product(s).**

There is no obligation to replace Products as the result of (i) accident, disaster or event of force majeure, (ii) misuse, fault or negligence of or by Buyer, (iii) use of the Products in a manner for which they were not designed, or (iv) improper storage and handling of the Products.

Current product instructions are available at [www.thermoscientific.com/pierce](http://www.thermoscientific.com/pierce). For a faxed copy, call 800-874-3723 or contact your local distributor.

© 2011 Thermo Fisher Scientific Inc. All rights reserved. Unless otherwise indicated, all trademarks are property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.