

## ACUITYAdvanced Biotin Free Polymer Kit (Previously Covance catalog# SIG-32906)

<b>Catalog# / Size</b>	931201 / 500 tests
<b>Regulatory Status</b>	RUO
<b>Other Names</b>	Detection Kit
<b>Previously</b>	Covance Catalog# SIG-32906

### Kit Contents

---

<b>Kit Contents</b>	<ul style="list-style-type: none"><li>• Reagent 1, Serum Block, 50 mL</li><li>• Reagent 2, Boost, 50 mL</li><li>• Reagent 3, HRP Polymer, 50 mL</li></ul>
---------------------	---

### Product Details

---

<b>Storage &amp; Handling</b>	Store between 2°C and 8°C.
<b>Application</b>	<a href="#">IHC</a>
<b>Recommended Usage</b>	Each lot of this product is quality control tested by immunohistochemistry.  Optimal working dilution and incubation times are to be determined by the investigator. See ACUITYAdvanced Protocol for more details.
<b>Application Notes</b>	ACUITYAdvanced is a biotin-free HRP polymer detection system for use with polyclonal and monoclonal antibodies made in rabbit and mouse. This technology provides excellent sensitivity and high specificity. ACUITYAdvanced is recommended for formalin fixed paraffin-embedded tissue sections. ACUITY Advanced is offered in a variety of kit sizes, and bulk volumes are available upon request.

### Antigen Details

---

<b>Gene ID</b>	NA
----------------	----

### Related Protocols

---

[AcuityADVANCED Protocol](#)

For research use only. Not for diagnostic use. Not for resale. BioLegend will not be held responsible for patent infringement or other violations that may occur with the use of our products.

\*These products may be covered by one or more Limited Use Label Licenses (see the BioLegend Catalog or our website, [www.biolegend.com/ordering#license](http://www.biolegend.com/ordering#license)). BioLegend products may not be transferred to third parties, resold, modified for resale, or used to manufacture commercial products, reverse engineer functionally similar materials, or to provide a service to third parties without written approval of BioLegend. By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.

BioLegend Inc., 8999 BioLegend Way, San Diego, CA 92121 [www.biolegend.com](http://www.biolegend.com)  
Toll-Free Phone: 1-877-Bio-Legend (246-5343) Phone: (858) 768-5800 Fax: (877) 455-9587