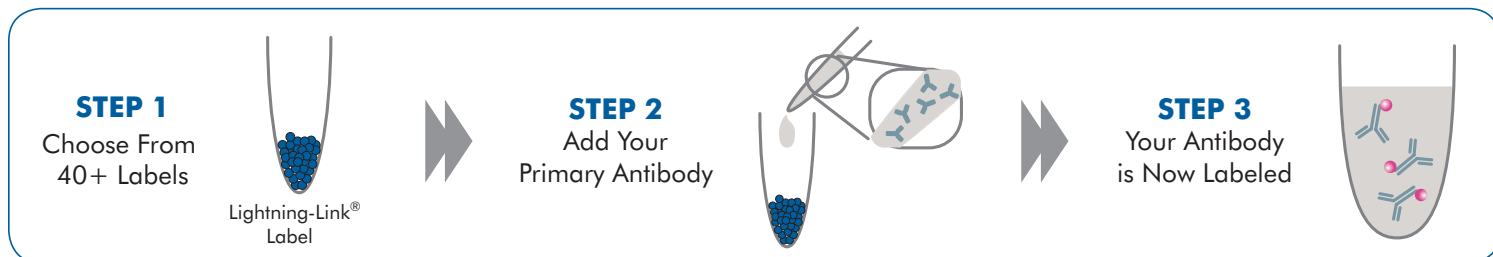


Novus Antibody Labeling Kits



- Label primary antibodies in 30 seconds
- Chose from 40 different labels
- Use with antibodies, proteins and peptides



Available Conjugates

Fluorescent Dyes

AMCA	FluorProbes® 594
Atto 390, 488, 565, 594, 633, 680, 700	FITC
Cyanine Dye Cy3, Cy5, Cy5.5	Rhodamine
DyLight™ 350, 405, 488, 550, 594, 633, 650, 680, 755, 800	Texas Red®

Fluorescent Proteins

Allophycocyanin (APC)	PerCP
B-Phycoerythrin (BPE)	R-Phycocyanin (RPC)

Tandem Dyes

Allophycocyanin/Cy5.5	PE/Cy5
Allophycocyanin/Cy7	PE/Cy5.5
PerCP/Cy5.5	PE/Cy7

Biotin/Streptavidin

Avidin	Biotin (Type A)
Streptavidin	Biotin (Type B)

Enzyme Conjugates

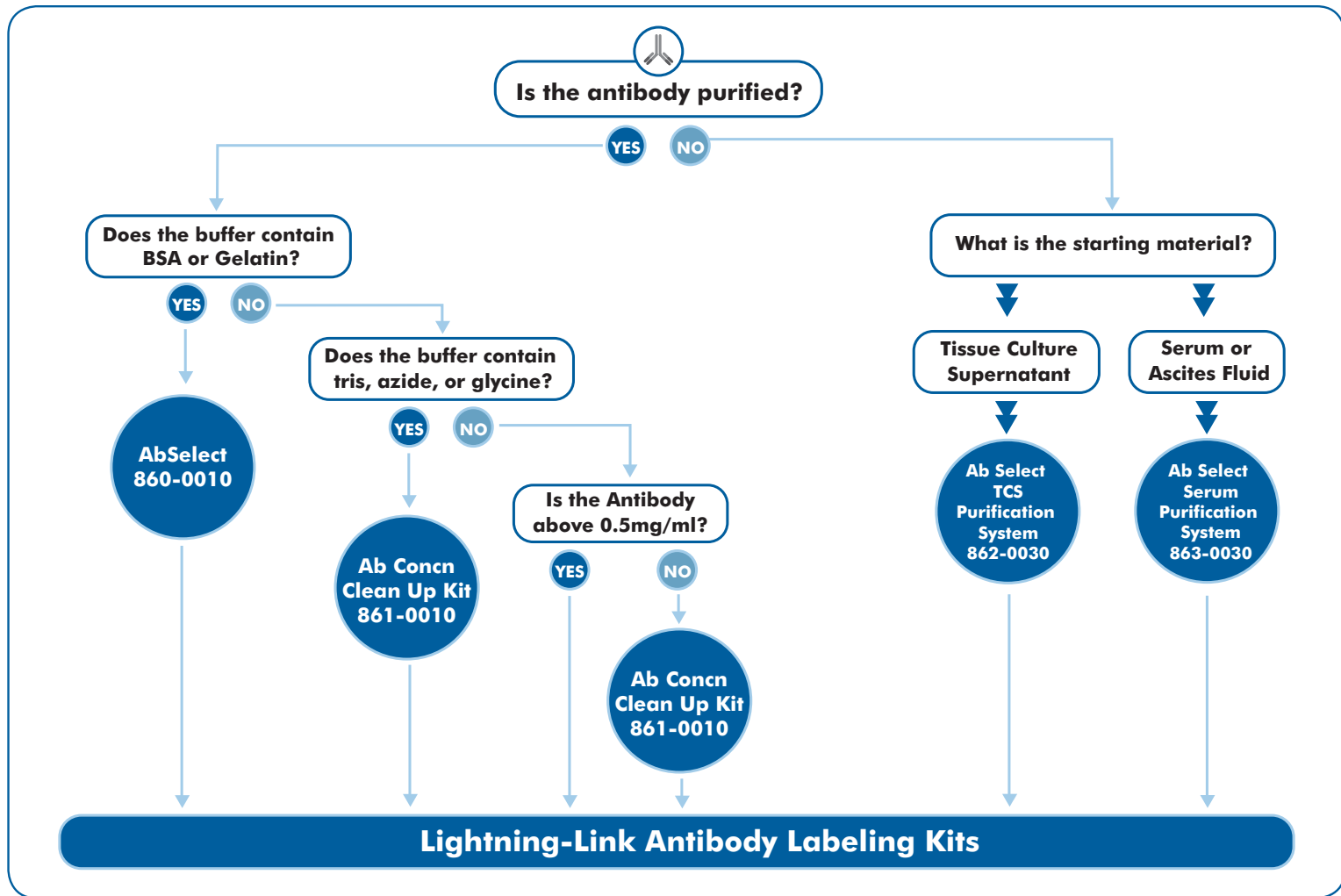
Alkaline Phosphatase (AP)
Glucose Oxidase (GOx)
Horseradish Peroxidase (HRP)

We also offer custom conjugation services. Our Antibody Lab will custom label many of the primary antibodies available on our website to over 18 different fluorescent or enzymatic labels.

For a complete product list, visit novusbio.com/labeling-kits

Lightning-Link® is a registered trademark of Innova Biosciences. DyLight™ is a trademark of Thermo Fisher Scientific Inc. Texas Red® is a registered trademark of Life Technologies Corporation. FluoroProbes® is a registered trademark of Interchim. ATTO Dyes are manufactured by ATTO-TEC GmbH. Cy5™ is a trademark of GE Healthcare.

Learn more | novusbio.com



AbSelect Antibody Purification System (860-0010)

Remove unwanted proteins, preservatives and amino acids that can interfere with labeling reactions. Purify antibodies from crude samples (ascites fluid or serum) and buffers containing high concentrations of BSA, azide or Tris. Purify up to 0.5 mg of antibody in a volume of 100 to 500 µl.

AbSelect Serum Purification Systems (863-0030)

Purify up to 20 mg of antibody from ascites fluid and serum.

AbSelect TCS System (862-0030)

Purify antibodies from hybridoma cell lines and supplied in tissue culture supernatant (TCS). Purify antibodies from 10-50 ml of TCS.

Antibody Concentration and Cleanup Kit (861-0010)

Allows for the quick and easy concentration of antibodies. The kit can also be used to reduce the concentration of any undesirable low molecular weight additives.

NOVUS USA

P: 303.730.1950
TF: 888.506.6887
F: 303.730.1966
E: novus@novusbio.com

R&D SYSTEMS EUROPE

P: +44 (0) 1235 529449
TF: 0800 37 34 15
F: +44 (0) 1235 533420
E: info.uk@bio-techne.com

NOVUS CANADA

P: 905.827.6400
TF: 855.668.8722
F: 905.827.6402
E: canada@novusbio.com

TECHNICAL SUPPORT

E: technical@novusbio.com

WORLDWIDE DISTRIBUTORS

www.novusbio.com/distributors

Novus Biologicals is a Bio-Techne brand.

Live Chat featured on our website, www.novusbio.com

Learn more | novusbio.com

bio-techne®

Global bio-techne.com info@bio-techne.com TEL +1 612 379 2956 North America TEL 800 343 7475
Europe | Middle East | Africa TEL +44 (0)1235 529449 China info.cn@bio-techne.com TEL +86 (21) 52380373

For research use or manufacturing purposes only. Trademarks and registered trademarks are the property of their respective owners.