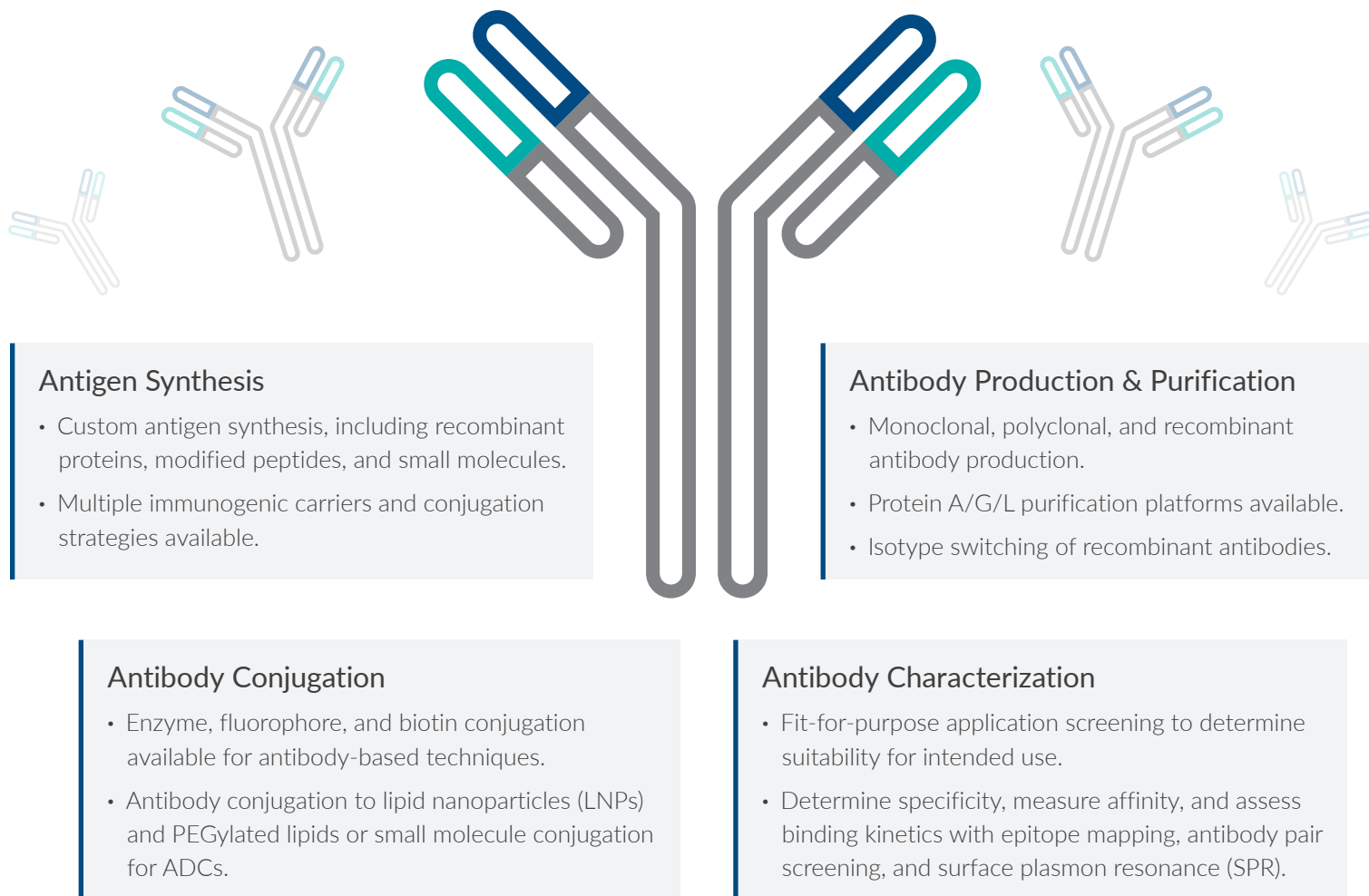


Antibody Services

Antibodies are essential for life science research because of their versatility for the detection, identification, and quantification of target molecules using a wide range of techniques. Cayman offers antibody services from antigen to application to support your research needs. Whether you need production of a novel antibody, synthesis of an antibody-drug conjugate (ADC), or assistance performing an antibody-based technique, Cayman has the expertise and facilities to advance your project to the next phase, step, or milestone.

Our Services for Antibody Production



Our Core Facilities Support Every Phase of Development

- Protein Production
- Cell & Molecular Biology
- LC-MS/MS
- Custom Assay Development
- Structural Biology
- Custom Synthesis

Our Services for Antibody-Based Applications

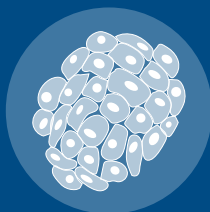
Our services support antibody-driven discoveries in applications ranging from target and complex isolation to expression-level analyses and effector antibodies in drug development.

Incorporate your custom antibody into your research or leverage our in-house expertise to perform these techniques for target isolation and sample analysis. We can also utilize your antibodies for fit-for-purpose application development/validation.



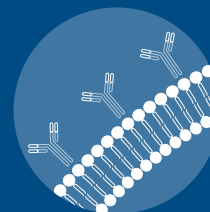
Target Isolation & Protein-Protein Interactions

- Immunoprecipitation from cells or tissue (antibody-coupled resin and sorbent development available)
- Mass spectrometry-based proteomics and immunopeptidome profiling



Expression-Level Analysis

- Western blot
- ELISA
- Flow cytometry
- Immunofluorescence
- Immunohistochemistry (IHC)



Effector Antibodies

- Conjugation of small molecule drugs for ADC development
- Addition to LNPs for targeting
- Isotype switching for recombinant production

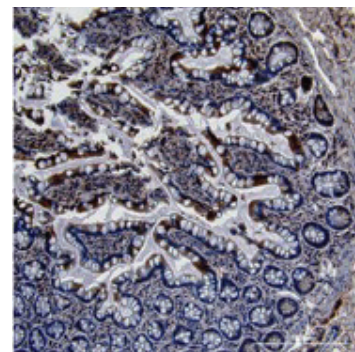
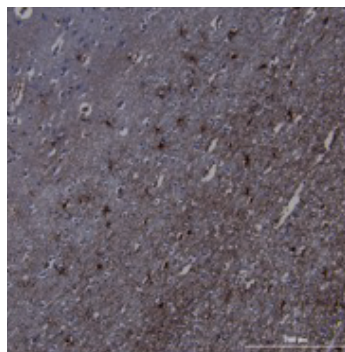
Scientific Poster

Development and Testing of IHC-Specific Antibodies at Cayman

Learn how our extensive experience with antigen and antibody development supports tissue sectioning and IHC testing on formalin-fixed, paraffin-embedded (FFPE) tissue blocks.

See the workflow at

www.caymanchem.com/IHC-poster



IHC imaging of GFAP in human cerebral cortex (left) and TMPRSS2 in human small intestine (right).

Learn more at www.caymanchem.com/antibody-services