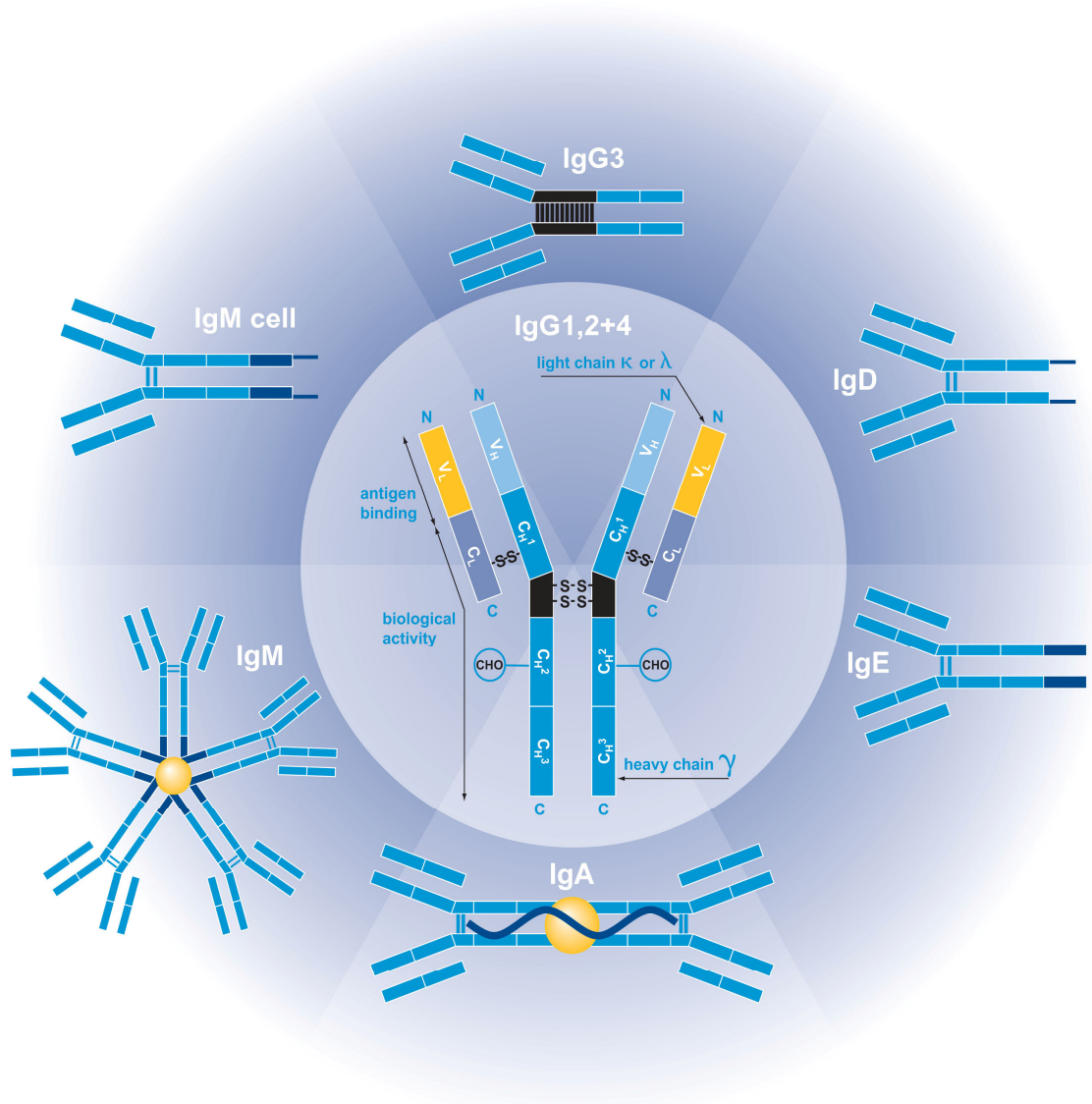


Antibody Isotypes



In mammals, antibodies can be divided into five isotypes or classes: IgG, IgM, IgA, IgD and IgE, based on the number of Y units and the type of heavy chain. Heavy chains of IgG, IgM, IgA, IgD, and IgE, are known as gamma, mu, alpha, delta, and epsilon, respectively.