

PNEUMOCOCCAL POLYSACCHARIDES – ADVANCING RESEARCH AND DEVELOPMENT

By Cara Wilder, PhD, ATCC, and Susan E. Witko, BS, Pfizer

Carbohydrates in the form of capsular polysaccharides and lipopolysaccharides constitute a major component of the cellular surface for many bacterial species.¹ In *Streptococcus pneumoniae*, capsular polysaccharides function as major virulence factors that protect the bacterium from phagocytosis by host immune cells. Within this species, capsular polysaccharides exhibit enormous structural diversity, resulting in significant differences in immunogenicity and antigenicity. In turn, this has presented numerous challenges toward the prevention and serological analysis of pneumococcal infections.² In the present article, we will discuss the importance of capsular polysaccharides in *S. pneumoniae* infection and will describe the use of purified serotype-specific polysaccharides in developing and evaluating pneumococcal vaccines, verifying novel immunoassays, and tracking bacterial disease epidemiology.

S. pneumoniae is a major human pathogen known to cause both invasive and non-invasive infections such as pneumonia, meningitis, bacteremia, otitis media, and sinusitis.^{2,3} More importantly, this bacterium is a leading cause of vaccine-preventable morbidity and mortality worldwide.⁴ In the United States, pneumococcal pneumonia is estimated to result in 150,000 hospitalizations annually with 5-7% of cases resulting in death.⁵ Moreover, the Global Disease Burden study from 2018 reported that *S. pneumoniae* was the leading cause of lower respiratory infection morbidity and mortality globally and contributed to more deaths than all other etiologies combined in 2016.⁶

This pneumococcal pathogen causes disease through the production of numerous virulence factors, including pneumolysin, pneumococcal surface proteins, and capsular polysaccharides.² Of these, the capsular polysaccharides—which are polymeric, surface-exposed carbohydrate molecules that encapsulate the bacteria—are the most important virulence factors as they shield the bacteria from neutrophil clearance. Currently, more than 95 different capsular serotypes have been identified, each distinguishable by serological response, variations in chemical structure, and related genetic mutations.^{2,7-13} These unique differences can associate with distinct epidemiological properties, including variations in carriage or prevalence of disease. For example, capsular types with higher ratios of charge to carbon may be physically larger, making them more resistant to neutrophil clearance.^{14,15}

Because of their importance in pneumococcal pathogenicity, capsular polysaccharides have been components in a number of serotype-specific vaccines. In 1983, a 23-valent pneumococcal polysaccharide vaccine (PPV23, Pneumovax23[®]; Merck Sharp & Dohme Corp.) was developed; this vaccine contained the pooled capsular polysaccharides purified from 23 different serotypes.¹⁶ Due to poor immunogenicity in infants, PPV23 was not approved for use in children younger than 2 years of age. In 2000, after nearly two decades without a suitable vaccine for juveniles, a 7-valent pneumococcal conjugate vaccine (7vPnC, Prevnar[®]; Wyeth Pharmaceuticals LLC, a subsidiary of Pfizer Inc.) was developed and licensed for use in infants and young children in the United States. This vaccine was created through the covalent coupling of a protein carrier to the purified polysaccharides from serotypes 4, 6B, 9V, 14, 18C, 19F, and 23F.¹⁷ To extend vaccine coverage to additional serotypes, a 13-valent pneumococcal conjugate vaccine (13vPnC, Prevnar 13[®]; Wyeth Pharmaceuticals LLC, a subsidiary of Pfizer Inc.) was later licensed for pediatric and adult use. 13vPnC is composed of the same serotype-conjugates included in 7vPnC as well as polysaccharide conjugates for serotypes 1, 3, 5, 6A, 7F, and 19A.^{18,19}

While widespread use of these vaccines has resulted in a reduction in pneumococcal pneumonia in targeted populations, non-vaccine serotypes have increased over time.²⁰ To address this problem, a 15-valent pneumococcal conjugate vaccine (PCV15, Vaxneuvance™; Merck Sharp & Dohme Corp.) and a 20-valent pneumococcal conjugate vaccine (20vPnC, Prevnar 20®; Wyeth Pharmaceuticals LLC, a subsidiary of Pfizer Inc.) were developed and recently approved for use in adults aged 18 years and older.²¹ These vaccines build upon the efficacy of 13vPnC through the incorporation of additional serotypes to provide further coverage for serotypes associated with pneumococcal disease worldwide. PCV15 includes serotypes 22F and 33F, which are among the serotypes known to cause invasive pneumococcal disease following the widespread use of 13vPnC.²¹ 20vPnC includes the addition of serotypes 8, 10A, 11A, 12F, 15B, 22F, and 33F, which have been associated with high case-fatality rates, antibiotic resistance, and meningitis.²²

In addition to their role in pneumococcal vaccine development, purified polysaccharides are also necessary for evaluating how effectively a vaccine can induce an immune response. Currently, the gold standard for measuring immunogenicity is via a standardized enzyme-linked immunosorbent assay (ELISA) protocol designed for the quantitation of human IgG antibodies specific for *S. pneumoniae* capsular polysaccharides.²³ The protocol for this assay, which was developed in 2000 by representatives from academia, government, industry, and the World Health Organization (WHO), is provided as a training manual containing the standard operating procedures for the preparation, execution, and analysis of the ELISA.²⁴ In this assay, ELISA plates are first coated with purified, serotype-specific capsular pneumococcal polysaccharides (ATCC, Manassas, VA). Following adsorption of the individual polysaccharides, dilutions of absorbed human sera are then added to the ELISA plates. The serotype-specific antibodies that remain bound to the purified pneumococcal polysaccharides are then detected and quantitated using an anti-human IgG antibody conjugated with alkaline phosphatase.²⁴

In recent years, capsular polysaccharides have been similarly used in the development and evaluation of novel immunoassays. Though the WHO ELISA protocol provides a standardized, accurate method for the evaluation of vaccine efficacy, it has come under scrutiny as it is laborious and consumes large volumes of serum. To this end, the development of novel multiplex immunoassays that allow for the simultaneous measurement of specific antibodies against various pneumococcal polysaccharides have been explored.²⁵⁻²⁹ For example, a 2012 study by Klein et al. described the development and characterization of a multiplex bead-based antibody quantification assay (MBIA) based on Luminex xMAP technology (Luminex, Austin, TX) that required the conjugation of purified serotype-specific capsular pneumococcal polysaccharides (ATCC, Manassas, VA) to fluorescent beads. In this assay, the polysaccharide-bead conjugate mixture was incubated with various dilutions of pre-adsorbed serum. Following incubation, bound antibodies were quantified using phycoerythrin-labeled goat anti-human reporter antibodies.²⁵ As compared to the traditional ELISA protocol, the MBIA procedure was able to provide a more efficient, cost-effective means of analyzing small volumes of serum.

Along with their function in detecting and enumerating anti-pneumococcal antibodies, the combinatorial use of capsular polysaccharides in immunoassays can be used for tracking the epidemiology of serotype-specific pneumococcal infections following vaccination. The gold-standard method for tracking invasive *S. pneumoniae* infection has occurred through traditional culture-based assays. However, good-quality samples are not always available, culturing and serological analyses can be fairly time consuming and laborious, and the detection of viable pneumococci can be hampered by use of antibiotics.^{30,31} To assess the positive impact of novel pneumococcal conjugate vaccines, serotype-specific diagnostics can provide valuable data on how effectively the vaccine reduces disease in the population and can assist in monitoring for the potential emergence of rare serotypes.

Pfizer Inc, a global leader in the biopharmaceutical industry committed to producing quality products that make an impact on health worldwide, is collaborating with ATCC, a leading provider of standards for the global scientific community, to ensure that researchers worldwide have access to purified, high-quality polysaccharides. The supply of pneumococcal polysaccharides is essential in continuing to drive pneumococcal research, predicting future epidemiological shifts, and combating this deadly pathogen.

Overall, capsular polysaccharides are bacterial antigens that play a significant role in pneumococcal pathogenicity. By targeting these polymeric carbohydrates using polysaccharide or polysaccharide conjugate vaccines, the disease burden of *S. pneumoniae* serotypes covered by these vaccines has been reduced significantly. Availability of purified pneumococcal polysaccharides has also facilitated the development of serological assays and diagnostics and improved the understanding of disease epidemiology.

REFERENCES

1. Weintraub A. Immunology of bacterial polysaccharide antigens. *Carbohydr Res* 338(23): 2539-2547, 2003. PubMed: [14670715](#)
2. Song JY, Nahm MH, Moseley MA. Clinical implications of pneumococcal serotypes: invasive disease potential, clinical presentations, and antibiotic resistance. *J Korean Med Sci* 28(1): 4-15, 2013. PubMed: [23341706](#)
3. Kaltoft MS, Skov Sorensen UB, Slotved HC, Konradsen HB. An easy method for detection of nasopharyngeal carriage of multiple *Streptococcus pneumoniae* serotypes. *J Microbiol Methods* 75(3): 540-544, 2008. PubMed: [18801391](#)
4. Frenkel LD. The global burden of vaccine-preventable infectious diseases in children less than 5 years of age: Implications for COVID-19 vaccination. How can we do better? *Allergy Asthma Proc* 42(5): 378-385, 2021. PubMed: [34474707](#)
5. World Health Organization. Pneumococcal Disease, <<https://www.nfd.org/infectious-diseases/pneumococcal/>>, 2022.
6. GBD 2016 Lower Respiratory Infections Collaborators. Estimates of the global, regional, and national morbidity, mortality, and aetiologies of lower respiratory infections in 195 countries, 1990–2016: a systematic analysis for the Global Burden of Disease Study 2016. *Lancet Infect Dis* 18(11): 1191-1210, 2018. PubMed: [30243584](#)
7. Park IH, et al. Discovery of a new capsular serotype (6C) within serogroup 6 of *Streptococcus pneumoniae*. *J Clin Microbiol* 45(4): 1225-1233, 2007. PubMed: [17267625](#)
8. Zartler ER, et al. Structure of the capsular polysaccharide of pneumococcal serotype 11A reveals a novel acetylglycerol that is the structural basis for 11A subtypes. *J Biol Chem* 284(11): 7318-7329, 2009. PubMed: [19114709](#)
9. Jin P, et al. First report of putative *Streptococcus pneumoniae* serotype 6D among nasopharyngeal isolates from Fijian children. *J Infect Dis* 200(9): 1375-1380, 2009. PubMed: [19803727](#)
10. Calix JJ, Dagan R, Pelton SI, Porat N, Nahm MH. Differential occurrence of *Streptococcus pneumoniae* serotype 11E between asymptomatic carriage and invasive pneumococcal disease isolates reflects a unique model of pathogen microevolution. *Clin Infect Dis* 54(6): 794-799, 2012. PubMed: [22267713](#)
11. Calix JJ, et al. Biochemical, genetic, and serological characterization of two capsule subtypes among *Streptococcus pneumoniae* Serotype 20 strains: discovery of a new pneumococcal serotype. *J Biol Chem* 287(33): 27885-27894, 2012. PubMed: [22736767](#)
12. Henrichsen J. Six newly recognized types of *Streptococcus pneumoniae*. *J Clin Microbiol* 33(10): 2759-2762, 1995. PubMed: [8567920](#)
13. Geno KA, et al. Pneumococcal Capsules and Their Types: Past, Present, and Future. *Clin Microbiol Rev* 28(3): 871-899, 2015. PubMed: [26085553](#)
14. Hausdorff WP, Feikin DR, Klugman KP. Epidemiological differences among pneumococcal serotypes. *Lancet Infect Dis* 5(2): 83-93, 2005. PubMed: [15680778](#)
15. Weinberger DM, et al. Pneumococcal capsular polysaccharide structure predicts serotype prevalence. *PLoS Pathog* 5(6): e1000476, 2009. PubMed: [19521509](#)
16. Robbins JB, et al. Considerations for formulating the second-generation pneumococcal capsular polysaccharide vaccine with emphasis on the cross-reactive types within groups. *J Infect Dis* 148(6): 1136-1159, 1983. PubMed: [6361173](#)
17. Advisory Committee on Immunization Practices. Preventing pneumococcal disease among infants and young children: recommendation of the Advisory Committee on Immunization Practices (ACIP). *MMWR Recomm Rep* 49(RR-9): 1-35, 2000. PubMed: [11055835](#)
18. US Food & Drug Administration. Product approval information---licensing action, package insert: Prevnar 13 (pneumococcal 13-valent conjugate vaccine), Pfizer. Rockville, MD: US Department of Health and Human Services, Food and Drug Administration., <<http://www.fda.gov/BiologicsBloodVaccines/Vaccines/ApprovedProducts/ucm201667.htm>>, 2010.
19. US Food & Drug Administration. Prevnar 13, <<https://www.fda.gov/vaccines-blood-biologics/vaccines/prevnar-13>>, 2019.
20. Stacey HL, et al. Safety and immunogenicity of 15-valent pneumococcal conjugate vaccine (PCV-15) compared to PCV-13 in healthy older adults. *Hum Vaccin Immunother* 15(3): 530-539, 2019. PubMed: [30648919](#)
21. Centers for Disease Control and Prevention. Use of 15-Valent Pneumococcal Conjugate Vaccine and 20-Valent Pneumococcal Conjugate Vaccine Among U.S. Adults: Updated Recommendations of the Advisory Committee on Immunization Practices — United States, 2022, <<https://www.cdc.gov/mmwr/volumes/71/wr/mm7104a1.htm>>, 2022.
22. Pfizer. U.S. FDA Approves PREVNAR 20™, Pfizer's Pneumococcal 20-valent Conjugate Vaccine for Adults Ages 18 Years or Older, <<https://www.pfizer.com/news/press-release/press-release-detail/us-fda-approves-prevnar-20tm-pfizers-pneumococcal-20-valent>>, 2021.
23. LaFon David C, Nahm MH. Measuring immune responses to pneumococcal vaccines. *J Immunol Methods* 461: 37-43, 2018. PubMed: [30098317](#)
24. World Health Organization Pneumococcal Serology Reference Laboratories at the Institute of Child Health, U. C. L., London, England, and the Department of Pathology at the University of Alabama at Birmingham, Birmingham Alabama, USA. Training manual for Enzyme linked immunosorbent assay for the quantification of *Streptococcus pneumoniae* serotype specific IgG (Pn PS ELISA). 007sp Version, 2000.

25. Klein DL, et al. Development and characterization of a multiplex bead-based immunoassay to quantify pneumococcal capsular polysaccharide-specific antibodies. *Clin Vaccine Immunol* 19(8): 1276-1282, 2012. PubMed: [22739691](#)
26. Goldblatt D, Ashton L, Zhang Y, Antonello J, Marchese RD. Comparison of a new multiplex binding assay versus the enzyme-linked immunosorbent assay for measurement of serotype-specific pneumococcal capsular polysaccharide IgG. *Clin Vaccine Immunol* 18(10): 1744-1751, 2011. PubMed: [21813660](#)
27. Elberse KE, Tcherniaeva I, Berbers GA, Schouls LM. Optimization and application of a multiplex bead-based assay to quantify serotype-specific IgG against *Streptococcus pneumoniae* polysaccharides: response to the booster vaccine after immunization with the pneumococcal 7-valent conjugate vaccine. *Clin Vaccine Immunol* 17(4): 674-682, 2010. PubMed: [20130129](#)
28. Pavliakova D, et al. Development and Validation of 13-plex Luminex-Based Assay for Measuring Human Serum Antibodies to *Streptococcus pneumoniae* Capsular Polysaccharides. *mSphere* 3(4): e00128-18, 2018. PubMed: [30089645](#)
29. Tan CY, et al. Evaluation of a Validated Luminex-Based Multiplex Immunoassay for Measuring Immunoglobulin G Antibodies in Serum to Pneumococcal Capsular Polysaccharides. *mSphere* 3(4): e00127-18, 2018. PubMed: [30089644](#)
30. Ewig S, et al. Factors associated with unknown aetiology in patients with community-acquired pneumonia. *Eur Resp J* 20(5): 1254-1262, 2002. PubMed: [12449182](#)
31. Musher DM, Montoya R, Wanahita A. Diagnostic value of microscopic examination of Gram-stained sputum and sputum cultures in patients with bacteremic pneumococcal pneumonia. *Clin Infect Dis* 39(2): 165-169, 2004. PubMed: [15307023](#)

Learn more about purified pneumococcal polysaccharides at
www.atcc.org/Polysaccharides

 10801 University Boulevard
Manassas, Virginia 20110-2209

 703.365.2700

 703.365.2701

 sales@atcc.org

 www.atcc.org

PP-072022-v05

©2022 American Type Culture Collection. The ATCC trademark and trade name, and any other trademarks listed in this publication are trademarks owned by the American Type Culture Collection unless indicated otherwise. *Streptococcus pneumoniae* bacteria image courtesy of CDC, Sarah Bailey Cutchin, and Dan Higgins.

These products are for laboratory use only. Not for human or diagnostic use. ATCC products may not be resold, modified for resale, used to provide commercial services or to manufacture commercial products without prior ATCC written approval.