



COLLEGE of AMERICAN  
PATHOLOGISTS

# CERTIFICATE OF ACCREDITATION

**Revvity Health Sciences Private  
Limited Laboratory  
Chennai, Tam, India  
Rajiv Rose, PhD**

CAP Number: 8649598  
AU-ID: 1653407

The organization named above meets all applicable standards for accreditation and is hereby accredited by the College of American Pathologists' Laboratory Accreditation Program. Reinspection should occur prior to **August 31, 2024** to maintain accreditation.

Accreditation does not automatically survive a change in director, ownership, or location and assumes that all interim requirements are met.

Michael Bradley Datto, MD, PhD, FCAP  
Chair, Accreditation Committee

Emily Volk, MD, FCAP  
President, College of American Pathologists

