

The College of American Pathologists
certifies that the laboratory named below

Asterand Inc
Biorepository
Detroit, Michigan
Chanel Smith, MS, BS

CAP Number: 8019367
AU-ID: 1628963

has met all applicable standards for accreditation and is hereby accredited by the
College of American Pathologists' Biorepository Accreditation Program. Reinspection
should occur prior to October 17, 2021 to maintain accreditation.

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



R. Bruce Williams, MD, FCAP

Chair, Accreditation Committee

President, College of American Pathologists