



COLLEGE of AMERICAN  
PATHOLOGISTS



The College of American Pathologists certifies  
that the laboratory named below

**Asterand Inc**  
**Biorepository**  
**Detroit, Michigan**  
**Charlenia Berry-Green, BA, AAS**

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, 2018 to maintain accreditation.

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

Chair, Commission on Laboratory Accreditation

President, College of American Pathologists