

Certificate of Registration

QUALITY MANAGEMENT SYSTEM - ISO 9001:2015

This is to certify that:

Asterand, Inc.
TechOne Building
440 Burroughs, Suite 501
Detroit
Michigan
48202
USA

Holds Certificate No:

FM 565646

and operates a Quality Management System which complies with the requirements of ISO 9001:2015 for the following scope:

The procurement and processing of high quality, well characterized human biospecimens for research use.

For and on behalf of BSI:



Carlos Pitanga, Chief Operating Officer Assurance – Americas

Original Registration Date: 2011-04-01

Latest Revision Date: 2018-08-24

Effective Date: 2017-04-01

Expiry Date: 2020-03-31

Page: 1 of 1



...making excellence a habit.™