

QUALITY MANAGEMENT SYSTEM CERTIFICATE

This certifies that the quality system of

BPS Bioscience, Inc.

6405 Mira Mesa Blvd., Unit #100, San Diego, CA 92121, USA

is registered by Systems Certification Body in recognition of a
Quality Management System which fulfills the requirement of

ISO 9001:2015

Scope of Registration

Design, expression, purification and characterization of active recombinant enzymes and
enzyme complexes for research applications

Certificate No: 1113189

Certificate Decision/Re-Issue Date:	02/14/2023
Certificate Issue Date:	09/28/2021
Certificate Expiry:	09/27/2024
Site Structure:	Single Site



SHIRLEY DEWI, SR. VICE PRESIDENT OF
MANAGEMENT SYSTEM REGISTRATION SERVICES

909.230.5526 | WWW.IAPMOSCB.ORG
5001 E. PHILADELPHIA ST, ONTARIO, CA 91761-2816

