

# Certificate of Registration



This is to certify that the Occupational Health & Safety Management System of:

**Cayman Pharma s.r.o.**

Práce 657, 277 11 Neratovice, Czech Republic

applicable to:

**Development and production of prostaglandins (active pharmaceutical ingredients) for human and veterinary use including associated activities**

has been assessed and registered by NQA against the provisions of:

**ISO 45001:2018**

This registration is subject to the company maintaining a quality management system, to the above standard, which will be monitored by NQA.

*NWright*

Managing Director



0015

Certificate No.	187218
Issue Date:	8 December 2019
Reissued:	5 December 2022
Valid Until:	7 December 2025
EAC Code:	13