

Certificate of Registration



This is to certify that the Environmental 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 14001:2015

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.	187217
Issue Date:	8 December 2007
Reissued:	5 December 2022
Valid Until:	7 December 2025
EAC Code:	13