

Separation of polycyclic aromatic hydrocarbons

HPTLC Silica gel 60 RP-18 F254s

Plate	HPTLC pre-coated plate RP-18 F 254s with concentrating zone (1.15498)
Eluent	Acetonitrile/methylene chloride/water 90/10/10 (v/v/v)
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Fluoranthene 2. Benzo(b)fluoranthene 3. Benzo(k)fluoranthene 4. Benzo(a)pyrene 5. Benzo(g,h, i)perylene 6. Indeno(1,2,3-cd)pyrene (all 0,05%)
Application volume	5 µl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 366 nm

