

Separation of some polynuclear aromatic hydrocarbons for analysis of water and environment

HPTLC Silica gel 60 RP-18 F254s

Plate	HPTLC pre-coated plate RP-18 F 254s with concentrating zone (1.15498)
Eluent	Acetonitrile/methanol 50/50 (v/v) temperature of development: -18°C
Migration Distance	6,5 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Benzo(e)pyrene 2. Chrysene 3. Triphenylene 4. Phenanthrene 5. Fluorene
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

