



Analysis of Sugars

CHROMATOGRAPHIC CONDITIONS

Column: Polypore CA (220 x 4.6 mm, 10 μ m) (07110095) with Polypore CA guard column (30 x 4.6 mm) (07110091)

Detection: RI

Mobile Phase: Water

Flow Rate: 0.2 mL/min. at 75° C, 100 psi

Sample: 10 μ L of standard or juice sample diluted 1:4 in water