

Product Data Sheet

Product Name: pVEC (Cadherin-5)

Catalog Number: AS-65465 (1 mg) Lot Number: See label on vial

Sequence: Leu-Leu-lle-lle-Leu-Arg-Arg-lle-Arg-Lys-Gln-Ala-His-Ala-his-Ser-Lys-OH

LLIILRRRIRKQAHAHSK

Molecular Weight: 2209.7

References:

Peptide Purity: Peak Area by HPLC ≥95%

Storage: This peptide is shipped at ambient temperature. Upon receipt, store

lyophilized peptide at -20°C or lower. Reconstituted peptide can be

aliquoted and stored at -20°C or lower.

Description: pVEC is a cell-penetrating peptide (CPP) that is derived from murine

vascular endothelian cadherin. It can permeate cell membrane at low micromolar concentration without significantly affecting cell membrane. pVEC can be conjugated with large molecules and used as a drug delivery

vehicle.

1. Nasrollahi, SA. et al. Chem Biol Drug Des **08**, 639-646 (2012).

2. Elmquist, A. et al. Exp Cell Res 269 (2), 237-244 (2001).

For Research Use Only