



## Product Data Sheet

---

<b>Product Name:</b>	Bovine $\beta$ -Casein, monophosphopeptide	
<b>Catalog Number:</b>	AS-61146 (1 mg)	Lot Number: See label on vial
<b>Sequence:</b>	H-Phe-Gln-pSer-Glu-Glu-Gln-Gln-Gln-Thr-Glu-Asp-Glu-Leu-Gln-Asp-Lys-OH (3-letter code) FQ-pS-EEQQQTEDELQDK (1-letter code)	
<b>Molecular Weight:</b>	2062.0	
<b>Peptide Purity:</b>	>95%	
<b>Appearance:</b>	Lyophilized white powder	

**Peptide Reconstitution:** Bovine  $\beta$ -Casein peptide is freely soluble in basic buffer.

**Storage:** Bovine  $\beta$ -Casein peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at  $-20^{\circ}\text{C}$  or lower. Reconstituted peptide can be aliquoted and stored at  $-20^{\circ}\text{C}$  or lower.

*For Research Use Only*