

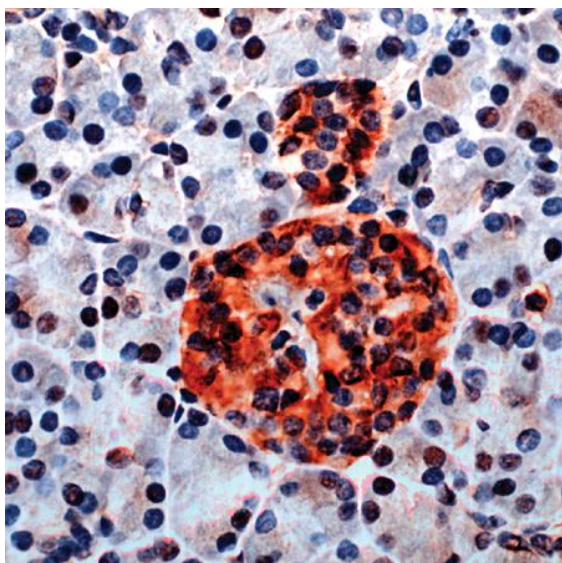


## Product Data Sheet

---

Product name:	<b>Anti-GAD-65 antibody, <i>polyclonal</i></b> ( <i>Glutamate decarboxylase 2</i> )	
Catalog number:	55990	
Lot number:	See label on vial	
Product description:	This antibody is supplied as 50 ug of affinity purified rabbit IgG in 1X PBS (pH 7.4) with 0.05% sodium azide.	
Immunogen:	Rabbit polyclonal anti-GAD65 antibody was raised against a synthetic peptide corresponding to the C-terminus of human GAD65 protein	
Species reactivity:	Confirmed: Human	
Validation:	IHC:	1: 50
	Other applications:	Not tested
Storage:	Store at 4°C for up to one year. For long term storage, aliquot and freeze. Avoid repeated freezing and thawing.	

This product is for *in vitro* research use only.



**Fig. 1.**  
IHC of human pancreas stained with  
polyclonal anti-GAD65 antibody  
(Cat # 55990)

**Target Information:**

Official symbol:	GAD2
Official full name:	Glutamate decarboxylase 2 (pancreatic islets and brain, 65kDa)
<i>Alternative Names:</i>	GAD65, 65 kDa glutamic acid decarboxylase, glutamate decarboxylase 65 kDa isoform
Organism:	Human
Length:	585 amino acids
Predicted MW:	65.4 kDa
Domains:	DDC / GAD / HDC / TyrDC pyridoxal-phosphate attachment site
Processing:	None identified
Subunit structure:	Homodimer
Tissue expression:	Not defined
Subcellular location:	Cytoplasm, cytosol, cytoplasmic vesicle, cell junction, synapse, presynaptic cell membrane, lipid-anchor, Golgi apparatus membrane, peripheral membrane. Note: Associated to cytoplasmic vesicles. In neurons, cytosolic leaflet of Golgi membranes and presynaptic clusters.
Function:	Gad65 catalyzes the production of GABA.
Disease association:	None identified

Source of target information is from HUGO Gene Nomenclature Committee, UniProt, ExPASy PROSITE.