

Murine Interferon Alpha Nomenclature

van Pesch <i>et al.</i> 2004			Hardy <i>et al.</i> 2004		Other Sequences	
Name	Accession	Activity	Name	Accession		
Alpha 1	AY225950 (C57Bl/6)	Mean	Alpha 1	NM_010502 (Balb/C)	AY226993 (129/Sv)	
Alpha 2	X01969 (Balb/C)	Mean	Alpha 2	NM_010503 (Balb/C)		
Alpha A	M28587 (Balb/C)	Mean	Alpha 3	M28587 (Balb/C)		
Alpha 4	X01973 (Balb/C)	High	Alpha 4	NM_010504 (Balb/C)	AY220463 (129/Sv)	
Alpha 5	X01971 (Balb/C)	Mean	Alpha 5	NM_010505 (Balb/C)	AY220464 (129/Sv)	
Alpha 6T	AY220465 (129/Sv)	Mean	Alpha 10		M23840 (Balb/C)	NM_026867 (129/Sv)
Alpha 7/10	M13710 (Swiss)	Low	Alpha 7	NM_088334 (C57Bl/6)	AY225952 (C57Bl/6)	
Alpha 8/6	X01972 (Balb/C)	Mean	Alpha 6	NM_008335	AY225953 (C57Bl/6)	NM_206871 (C57Bl/6)
Alpha 9	M13660 (Balb/C)	Mean	Alpha 9	M13660 (Balb/C)		
Alpha 11	M68944 (Swiss)	High	Alpha 11	NM_008333 (Swiss)	AY225954 (C57Bl/6)	
Alpha 12	AY225951 (C57Bl/6)	Mean	Alpha 12	AY190046 (C57Bl/6)		
Alpha 13	AY220461 (129/Sv)	Mean	Alpha 13	AY190047 (C57Bl/6)	NM_177347 (C57Bl/6)	
Alpha 14	AY220462 (129/Sv)	Mean				
Alpha B	L38698 (Balb/C)	Mean	Alpha 8	NM_008336		

Two groups have recently attempted to devise a unified nomenclature of the mouse IFN alpha subtypes. These two systems are very similar but with a few differences. We have chosen to use the van Pesch et al. nomenclature since this group expressed all the subtypes and attempted to determine relative antiviral activity, and they identified a new subtype, Alpha 14.