

## Human Interferon-Alpha Nomenclature

Product #	Nomenclature		Gene	Molecular Weight	Theoretical pI
11100-1	$\alpha$ A	$\alpha$ 2a	IFNA2	19241	5.99
11105-1	$\alpha$ 2	$\alpha$ 2b	IFNA2	19269	5.99
11115-1	$\alpha$ B2	$\alpha$ 8	IFNA8	19484	5.18
11120-1	$\alpha$ C	$\alpha$ 10	IFNA10	19406	5.70
11125-1	$\alpha$ D [Val 114]	$\alpha$ 1	IFNA1	19414	5.18
11130-1	$\alpha$ F	$\alpha$ 21	IFNA21	19312	5.99
11135-1	$\alpha$ G	$\alpha$ 5	IFNA5	19695	5.46
11145-1	$\alpha$ H2	$\alpha$ 14	IFNA14	19708	6.38
11150-1	$\alpha$ I	$\alpha$ 17	IFNA17	19299	5.46
11160-1	$\alpha$ J1	$\alpha$ 7	IFNA7	19606	5.87
11165-1	$\alpha$ K	$\alpha$ 6	IFNA6	20061	6.43
11175-1	$\alpha$ 1 [Ala114]	$\alpha$ D	IFNA1	19386	5.18
11177-1	$\alpha$ 4a	$\alpha$ M1	IFNA4	19379	5.58
11180-1	$\alpha$ 4b	$\alpha$ 4	IFNA4	19379	5.76
11190-1	$\alpha$ WA	$\alpha$ 16	IFNA16	19282	5.96

For more information of the various alleles of IFN alpha 2a, 2b, 4a, 4b, 1 and D

**References:**

1. Gewart, DR., et al. (1995). *J Interferon Cytokine Res.* 15 (5): 403
2. Hussain, M., et al. (1997). *J Interferon Cytokine Res.* 17 (9): 559
3. Hussian, M., et al. (2000). *J Interferon Cytokine Res.* 20: 763